

Covance Price List				
CatalogNo	Product Name	Size	EUR	Product Line
A488-101L-0100	HA.11 Monoclonal Antibody, Alexa Fluor® Labeled	0.1 mL	479	Covance Core
A488-114L-0100	Giantin Polyclonal Antibody, Alexa Fluor® Labeled	0.1 mL	479	Covance Core
A488-120L-0100	Alexa Fluor Labeled MAb414	0.1 ml	479	Added April 10th
A488-128L-0100	RNA Polymerase II CTD4H8 Monoclonal Antibody, Alexa Fluor® Labeled	0.1 mL	479	Covance Core
A488-150L-0100	9E10 Monoclonal Antibody (anti-c-myc), Alexa Fluor® Labeled	0.1 mL	479	Covance Core
A488-164L-0100	9E11 (anti c-myc) Monoclonal Antibody Alexa Fluor® Labeled	0.1 mL	479	Covance Core
A488-435L-0100	Neuronal Class III & beta;-Tubulin (TUJ1) Monoclonal Antibody, Alexa Fluor® Labeled	0.1 mL	479	Covance Core
A594-101L-0100	HA.11 Monoclonal Antibody, Alexa Fluor® Labeled	0.1 mL	479	Covance Core
AFC-101P-1000	HA.11 Monoclonal Antibody, Affinity Matrix	1.0 mL	667	Covance Core
AFC-115P-1000	GLU-GLU Monoclonal Antibody, Affinity Matrix	1.0 mL	667	Covance Core
AFC-130P-1000	AU1 Monoclonal Antibody, Affinity Matrix	1.0 mL	667	Covance Core
AFC-135P-1000	AU5 Monoclonal Antibody, Affinity Matrix	1.0 mL	667	Covance Core
AFC-150P-1000	9E10 Monoclonal Antibody, Affinity Matrix	1.0 mL	667	Covance Core
AFC-156P-1000	6-His Monoclonal Antibody, Affinity Matrix	1.0 mL	667	Covance Core
BIOT-101L-0100	HA.11 Monoclonal Antibody, Biotin Labeled	0.1 mL	467	Covance Core
BIOT-106L-0100	HA.11 Monoclonal Antibody, Biotin Labeled	0.1 mL	467	Covance Core
BIOT-115L-0100	GLU-GLU Monoclonal Antibody, Biotin Labeled	0.1 mL	467	Covance Core
BIOT-130L-0100	AU1 Monoclonal Antibody, Biotin Labeled	0.1 mL	467	Covance Core
BIOT-135L-0100	AU5 Monoclonal Antibody, Biotin Labeled	0.1 mL	467	Covance Core
BIOT-150L-0100	9E10 Monoclonal Antibody, Biotin Labeled	0.1 mL	467	Covance Core
BIOT-156L-0100	6-His Monoclonal Antibody, Biotin Labeled	0.1 mL	467	Covance Core
FITC-101L-0100	HA.11 Monoclonal Antibody, FITC Labeled	0.1 mL	467	Covance Core
FITC-106L-0100	Cre Recombinase Monoclonal Antibody, FITC Labeled	0.1 mL	467	Covance Core
FITC-115L-0100	GLU-GLU Monoclonal Antibody, FITC Labeled	0.1 mL	467	Covance Core
FITC-130L-0100	AU1 Monoclonal Antibody, FITC Labeled	0.1 mL	467	Covance Core
FITC-135L-0100	AU5 Monoclonal Antibody, FITC Labeled	0.1 mL	467	Covance Core
FITC-150L-0100	9E10 Monoclonal Antibody, FITC Labeled	0.1 mL	467	Covance Core
FITC-156L-0100	6-His Monoclonal Antibody, FITC Labeled	0.1 mL	467	Covance Core
HRP-164L-0100	9E11 (anti c-myc) Monoclonal Antibody, HRP Labeled	0.1 mL	467	Covance Core
MMS-101P-1000	HA.11 Monoclonal Antibody, Purified	1.0 mL	882	Covance Core
MMS-101P-0200	HA.11 Monoclonal Antibody, Purified	0.2 mL	400	Covance Core
MMS-101P-0500	HA.11 Monoclonal Antibody, Purified	0.5 mL	600	Covance Core
MMS-101R-1000	HA.11 Monoclonal Antibody	1.0 mL	773	Covance Core
MMS-101R-0200	HA.11 Monoclonal Antibody	0.2 mL	326	Covance Core
MMS-101R-0500	HA.11 Monoclonal Antibody	0.5 mL	498	Covance Core
MMS-102P-0100	NUP 153 Nuclear Pore Complex Protein Monoclonal Antibody	0.1 mL	348	Covance Core
MMS-102P-0500	NUP 153 Nuclear Pore Complex Protein Monoclonal Antibody	0.5 mL	655	Covance Core
MMS-106P-0200	Cre Recombinase Monoclonal Antibody, Purified	0.2 mL	400	Covance Core
MMS-106R-0200	Cre Recombinase Monoclonal Antibody	0.2 mL	326	Covance Core
MMS-107P-0100	Lamin A & C (XB10) Monoclonal Antibody, Purified	0.1 mL	348	Covance Core
MMS-110R-0200	Mannosidase II Monoclonal Antibody	0.2 mL	304	Covance Core
MMS-110R-0500	Mannosidase II Monoclonal Antibody	0.5 mL	482	Covance Core
MMS-112P-0200	GST Monoclonal Antibody, Purified	0.2 mL	400	Covance Core
MMS-112P-0500	GST Monoclonal Antibody, Purified	0.5 mL	600	Covance Core
MMS-112R-0200	GST Monoclonal Antibody	0.2 mL	326	Covance Core
MMS-112R-0500	GST Monoclonal Antibody	0.5 mL	498	Covance Core
MMS-115P-0200	GLU-GLU Monoclonal Antibody, Purified	0.2 mL	400	Covance Core
MMS-115P-0500	GLU-GLU Monoclonal Antibody, Purified	0.5 mL	600	Covance Core
MMS-115R-0200	GLU-GLU Monoclonal Antibody	0.2 mL	326	Covance Core
MMS-115R-0500	GLU-GLU Monoclonal Antibody	0.5 mL	498	Covance Core
MMS-116P-0250	HIV TAT Monoclonal Antibody, Purified	0.25 mL	422	Covance Core
MMS-118P-0200	Green Fluorescent Protein (GFP) Monoclonal Antibody, Purified	0.2 mL	400	Covance Core
MMS-118P-0500	Green Fluorescent Protein (GFP) Monoclonal Antibody, Purified	0.5 mL	600	Covance Core
MMS-118R-0200	Green Fluorescent Protein (GFP) Monoclonal Antibody	0.2 mL	326	Covance Core
MMS-118R-0500	Green Fluorescent Protein (GFP) Monoclonal Antibody	0.5 mL	498	Covance Core
MMS-120P-0100	Nuclear Pore Complex Proteins (MAb414) Monoclonal Antibody, Purified	0.1 mL	339	Covance Core
MMS-120P-0500	Nuclear Pore Complex Proteins (MAb414) Monoclonal Antibody, Purified	0.5 mL	571	Covance Core
MMS-120R-0200	Nuclear Pore Complex Proteins (MAb414) Monoclonal Antibody	0.2 mL	386	Covance Core
MMS-120R-0500	Nuclear Pore Complex Proteins (MAb414) Monoclonal Antibody	0.5 mL	482	Covance Core
MMS-125P-0200	KT3 Monoclonal Antibody, Purified	0.2 mL	400	Covance Core
MMS-125P-0500	KT3 Monoclonal Antibody, Purified	0.5 mL	600	Covance Core
MMS-125R-0200	KT3 Monoclonal Antibody	0.2 mL	326	Covance Core
MMS-125R-0500	KT3 Monoclonal Antibody	0.5 mL	498	Covance Core
MMS-126R-0200	RNA Polymerase II 8WG16 Monoclonal Antibody	0.2 mL	374	Covance Core
MMS-126R-0500	RNA Polymerase II 8WG16 Monoclonal Antibody	0.5 mL	498	Covance Core
MMS-128P-0100	RNA Polymerase II CTD4H8 Monoclonal Antibody	0.1 mL	400	Covance Core
MMS-129R-0200	RNA Polymerase II H5 Monoclonal Antibody	0.2 mL	374	Covance Core
MMS-129R-0500	RNA Polymerase II H5 Monoclonal Antibody	0.5 mL	498	Covance Core
MMS-130P-0200	AU1 Monoclonal Antibody, Purified	0.2 mL	400	Covance Core
MMS-130P-0500	AU1 Monoclonal Antibody, Purified	0.5 mL	600	Covance Core
MMS-130R-0200	AU1 Monoclonal Antibody	0.2 mL	326	Covance Core
MMS-130R-0500	AU1 Monoclonal Antibody	0.5 mL	498	Covance Core
MMS-134R-0200	RNA Polymerase II H14 Monoclonal Antibody	0.2 mL	374	Covance Core
MMS-134R-0500	RNA Polymerase II H14 Monoclonal Antibody	0.5 mL	498	Covance Core
MMS-135P-0200	AU5 Monoclonal Antibody, Purified	0.2 mL	400	Covance Core
MMS-135P-0500	AU5 Monoclonal Antibody, Purified	0.5 mL	600	Covance Core

CatalogNo	Product Name	Size	EUR	Product Line
MMS-135R-0200	AU5 Monoclonal Antibody	0.2 mL	326	Covance Core
MMS-135R-0500	AU5 Monoclonal Antibody	0.5 mL	498	Covance Core
MMS-137P-0250	Cytokeratin (AE1/AE3) Monoclonal Antibody	0.25 mL	281	Covance Core
MMS-139S-0250	Bromodeoxyuridine (BrdU) Monoclonal Antibody	0.25 mL	361	Covance Core
MMS-141P-0250	Rat VCAM-1 Monoclonal Antibody, Purified	0.25 mL	422	Covance Core
MMS-143R-0100	Ezrin Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-146P-0250	TNF-α Monoclonal Antibody, Purified	0.25 mL	400	Covance Core
MMS-147P-0100	GM130 Monoclonal Antibody	0.1 mL	400	Covance Core
MMS-148S-0250	Cytokeratin 7 (OV-TL 12/30) Monoclonal Antibody	0.25 mL	315	Covance Core
MMS-150P-1000	9E10 Monoclonal Antibody (anti-c-myc), Purified	1.0 mL	882	Covance Core
MMS-150P-0200	9E10 Monoclonal Antibody (anti-c-myc), Purified	0.2 mL	400	Covance Core
MMS-150P-0500	9E10 Monoclonal Antibody (anti-c-myc), Purified	0.5 mL	600	Covance Core
MMS-150R-1000	9E10 Monoclonal Antibody (anti-c-myc)	1.0 mL	773	Covance Core
MMS-150R-0200	9E10 Monoclonal Antibody (anti-c-myc)	0.2 mL	326	Covance Core
MMS-150R-0500	9E10 Monoclonal Antibody (anti-c-myc)	0.5 mL	498	Covance Core
MMS-151S-0250	Cytokeratin (MNF116) Monoclonal Antibody	0.25 mL	308	Covance Core
MMS-156P-0200	6-His Monoclonal Antibody, Purified	0.2 mL	400	Covance Core
MMS-156P-0500	6-His Monoclonal Antibody, Purified	0.5 mL	600	Covance Core
MMS-156R-0200	6-His Monoclonal Antibody	0.2 mL	326	Covance Core
MMS-156R-0500	6-His Monoclonal Antibody	0.5 mL	498	Covance Core
MMS-157S-0250	Cytokeratin 10/13 (DE-K13) Monoclonal Antibody	0.25 mL	330	Covance Core
MMS-158S-0250	Cytokeratin 19 (RCK 108) Monoclonal Antibody	0.25 mL	315	Covance Core
MMS-159S-0250	Cytokeratin 10 (DE-K10) Monoclonal Antibody	0.25 mL	417	Covance Core
MMS-161S-0250	Cytokeratin 17 (E3) Monoclonal Antibody	0.25 mL	404	Covance Core
MMS-162P-0250	Human Cytokeratin-8 (HK-8) Monoclonal Antibody, Purified	0.25 mL	489	Covance Core
MMS-164P-0100	9E11 (anti c-myc) Monoclonal Antibody, Purified	0.1 mL	400	Covance Core
MMS-168P-0100	Beta-Galactosidase Monoclonal Antibody	0.1 mL	400	Covance Core
MMS-172R-0500	L1 NCAM Monoclonal Antibody	0.5 mL	482	Covance Core
MMS-174P-0250	Cytokeratin 20 (Ks 20.8) Monoclonal Antibody	0.25 mL	368	Covance Core
MMS-176P-0100	Thrombomodulin (Phx-01) Monoclonal Antibody, Purified	0.1 mL	400	Covance Core
MMS-178P-0100	Thrombomodulin (Phx-02) Monoclonal Antibody, Purified	0.1 mL	400	Covance Core
MMS-181P-0250	Cytokeratin 5/6 (D5/16 B4) Monoclonal Antibody	0.25 mL	650	Covance Core
MMS-187P-0250	HIV gp41 Monoclonal Antibody, Purified	0.25 mL	422	Covance Core
MMS-188P-0200	Kinesin Heavy Chain Monoclonal Antibody, Purified	0.2 mL	400	Covance Core
MMS-193P-0250	HIV gp120 Monoclonal Antibody, Purified	0.25 mL	422	Covance Core
MMS-196R-0250	SR Proteins and Splicing Factors, Monoclonal Antibody	0.25 mL	400	Covance Core
MMS-197R-0250	SR Proteins and Splicing Factors, Monoclonal Antibody	0.25 mL	400	Covance Core
MMS-198P-0200	Kinesin II Monoclonal Antibody	0.2 mL	400	Covance Core
MMS-201R-0100	HOX C4 Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-228R-0100	TATA Binding Proteins (TBP) Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-233R-0100	TATA Binding Proteins (TBP) Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-248R-0100	O-GlcNAc Monoclonal Antibody	0.1 mL	298	Covance Core
MMS-248R-0500	O-GlcNAc Monoclonal Antibody	0.5 mL	498	Covance Core
MMS-257P-0200	Ubiquitin (P4D1) Monoclonal Antibody	0.2 mL	498	Covance Core
MMS-258R-0200	Ubiquitin (P4G7) Monoclonal Antibody	0.2 mL	326	Covance Core
MMS-258R-0500	Ubiquitin (P4G7) Monoclonal Antibody	0.5 mL	498	Covance Core
MMS-259R-0100	E1 Ubiquitin Activating Enzyme, Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-261R-0100	Msx-1 Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-261R-0500	Msx-1 Monoclonal Antibody	0.5 mL	498	Covance Core
MMS-263R-0100	Ku, Small Subunit Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-264R-0100	Ubiquitin Ubi-1 Monoclonal Antibody	0.1 mL	356	Covance Core
MMS-267R-0100	Ku, Large Subunit Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-269R-0100	Chromatin Assembly Factor 1 (CAF-1) Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-273R-0100	Transcription Factor IIB (TFIIB) Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-282R-0100	Insulin Degrading Enzyme (IDE) Monoclonal Antibody	0.1 mL	298	Covance Core
MMS-286R-0100	HOX C8 [C952-12E] Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-286R-0500	HOX C8 [C952-12E] Monoclonal Antibody	0.5 mL	467	Covance Core
MMS-318P-0100	Sodium Calcium Exchanger 1 Monoclonal Antibody	0.1 mg	429	Covance Core
MMS-359S-1000	Presenilin 2 Monoclonal Antibody	1.0 mL	374	Covance Core
MMS-400P-0100	Cytoplasmic Dynein Monoclonal Antibody, Purified	0.1 mL	297	Covance Core
MMS-400R-0100	Cytoplasmic Dynein Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-403P-0250	Phosphotyrosine (PY20) Monoclonal Antibody, Purified	0.25 mL	400	Covance Core
MMS-405P-0250	Phosphotyrosine (6G9) Monoclonal Antibody, Purified	0.25 mL	400	Covance Core
MMS-407R-0200	α-Tubulin Monoclonal Antibody	0.2 mL	422	Covance Core
MMS-408P-0250	Phosphotyrosine (1G2) Monoclonal Antibody, Purified	0.25 mL	400	Covance Core
MMS-410P-0250	β-Tubulin Monoclonal Antibody, Purified	0.25 mL	438	Covance Core
MMS-411R-0200	β-Tubulin Monoclonal Antibody	0.2 mL	422	Covance Core
MMS-412P-0250	Phosphotyrosine (2G8) Monoclonal Antibody, Purified	0.25 mL	400	Covance Core
MMS-413R-0200	Acetylated α-Tubulin, Monoclonal Antibody	0.2 mL	422	Covance Core
MMS-414P-0250	Phosphotyrosine (PY72) Monoclonal Antibody, Purified	0.25 mL	400	Covance Core
MMS-416R-0100	Cardiac Troponin-I Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-418R-0100	Cardiac Troponin-I [Ser42] Monoclonal Antibody, Phospho-Specific	0.1 mL	297	Covance Core
MMS-422P-0200	Class II β-Tubulin (7B9) Monoclonal Antibody, Purified	0.2 mL	400	Covance Core
MMS-423P-0100	Clathrin Light Chain Monoclonal Antibody, Purified	0.1 mL	348	Covance Core
MMS-424P-0200	Class II β-Tubulin (6B1) Monoclonal Antibody, Purified	0.2 mL	400	Covance Core
MMS-427P-0100	Clathrin Heavy Chain Monoclonal Antibody, Purified	0.1 mL	348	Covance Core
MMS-429P-0100	Human Flightless-I Protein Monoclonal Antibody, Purified	0.1 mL	372	Covance Core

CatalogNo	Product Name	Size	EUR	Product Line
MMS-429R-0200	Human Flightless-I Protein Monoclonal Antibody	0.2 mL	400	Covance Core
MMS-434R-0100	Kaptein (Actin-Associated Protein 2E4) Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-435P-0100	Neuronal Class III β -Tubulin (TUJ1) Monoclonal Antibody, Purified	0.1 mL	361	Covance Core
MMS-435P-0250	Neuronal Class III β -Tubulin (TUJ1) Monoclonal Antibody, Purified	0.25 mL	438	Covance Core
MMS-438I-6000	Melanoma (HMB45) Monoclonal Antibody (Ready to Use)	6.0 mL	361	Covance Core
MMS-438P-0500	Melanoma (HMB45) Monoclonal Antibody	0.5 mL	436	Covance Core
MMS-442P-0100	Nonmuscle Myosin II Heavy Chain Monoclonal Antibody, Purified	0.1 mL	429	Covance Core
MMS-448P-0100	mAKAP (OR017.120) Monoclonal Antibody	0.1 mL	429	Covance Core
MMS-454S-0250	Desmin (D33) Monoclonal Antibody	0.25 mL	299	Covance Core
MMS-456S-0100	Myosin II (Pan) Monoclonal Antibody	0.1 mL	386	Covance Core
MMS-459S-0250	Vimentin (V9) Monoclonal Antibody	0.25 mL	288	Covance Core
MMS-460R-0100	Myosin Heavy Chain Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-461P-0500	Smooth Muscle Actin (CGA7) Monoclonal Antibody, Purified	0.5 mL	436	Covance Core
MMS-464S-0250	Vimentin (Vim 3B4) Monoclonal Antibody	0.25 mL	375	Covance Core
MMS-466S-0250	Smooth Muscle Actin (1A4) Monoclonal Antibody	0.25 mL	397	Covance Core
MMS-467S-0250	Sarcomeric Actin (Alpha Sr-1) Monoclonal Antibody	0.25 mL	376	Covance Core
MMS-468P-0500	Muscle Actin (HHF35) Monoclonal Antibody	0.5 mL	436	Covance Core
MMS-470R-0100	CD146 Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-470R-0500	CD146 Monoclonal Antibody	0.5 mL	535	Covance Core
MMS-472R-0100	Integrin β 1 Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-474R-0100	Integrin α 5 β 5 Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-474R-0500	Integrin α 5 β 5 Monoclonal Antibody	0.5 mL	535	Covance Core
MMS-476R-0100	VCAM-1 Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-480R-0100	Integrin β 1 (Activating) Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-482R-0100	E-Selectin Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-484R-0100	PECAM Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-484R-0500	PECAM Monoclonal Antibody	0.5 mL	535	Covance Core
MMS-485P-0100	MAP2 Monoclonal Antibody, Purified	0.1 mL	372	Covance Core
MMS-485R-0200	MAP2 Monoclonal Antibody	0.2 mL	386	Covance Core
MMS-486R-0100	P-Selectin Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-487P-0100	Detyrosinated α -Tubulin Monoclonal Antibody, Purified	0.1 mL	372	Covance Core
MMS-487R-0200	Detyrosinated α -Tubulin Monoclonal Antibody	0.2 mL	400	Covance Core
MMS-488R-0100	Integrin α 2 β 1 Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-489P-0100	α -Tubulin Monoclonal Antibody, Purified	0.1 mL	372	Covance Core
MMS-489R-0200	α -Tubulin Monoclonal Antibody	0.2 mL	400	Covance Core
MMS-490R-0100	Integrin α 3 β 1 Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-492R-0100	Integrin α 4 β 1 Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-493R-0100	MAP 2 Monoclonal Antibody	0.1 mL	360	Covance Core
MMS-494R-0100	Integrin β 2 Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-495R-0100	MAP 1B Monoclonal Antibody	0.1 mL	360	Covance Core
MMS-496R-0100	Integrin α 5 β 1 Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-0500P-0500	GFAP (GF2) Monoclonal Antibody	0.5 mL	436	Covance Core
MMS-501P-015	NMDA Receptor NR1, Monoclonal Antibody, Purified	0.015 mg	416	Covance Core
MMS-503P-0100	Tyrosine Hydroxylase Monoclonal Antibody	0.1 mL	436	Covance Core
MMS-518P-0200	Neuron-Specific Enolase (NSE-P1) Monoclonal Antibody	0.2 mL	400	Covance Core
MMS-519P-0200	Neuron-specific Enolase (NSE-P2) Monoclonal Antibody	0.2 mL	400	Covance Core
MMS-520R-0500	Tau 13 Monoclonal Antibody	0.5 mL	467	Covance Core
MMS-529R-0200	Synuclein- α / β Monoclonal Antibody	0.2 mL	422	Covance Core
MMS-530R-0200	Synuclein- α , Monoclonal Antibody	0.2 mL	422	Covance Core
MMS-531R-0200	Tau 46 Monoclonal Antibody	0.2 mL	422	Covance Core
MMS-532R-0200	Amyloid- β ; (A β) Monoclonal Antibody	0.2 mL	498	Covance Core
MMS-544R-0200	Neurofilament 160/200 kD (NF-M+H) Monoclonal Antibody	0.2 mL	422	Covance Core
MMS-545R-0200	Tau [Thr231] Monoclonal Antibody, Phospho-Specific	0.2 mL	422	Covance Core
MMS-546R-0200	Tau [Ser396] Monoclonal Antibody, Phospho-Specific	0.2 mL	422	Covance Core
MMS-555R-0200	Synuclein- α / β ; Nitrated Monoclonal Antibody	0.2 mL	422	Covance Core
MMS-556R-0200	Synuclein- α , Nitrated Monoclonal Antibody	0.2 mL	422	Covance Core
MMS-557R-0200	14-3-3 Monoclonal Antibody	0.2 mL	422	Covance Core
MMS-561R-0200	DJ-1 Monoclonal Antibody	0.2 mL	422	Covance Core
MMS-565R-0100	Bax Universal Monoclonal Antibody	0.1 mL	407	Covance Core
MMS-566R-0100	Bcl-XL Monoclonal Antibody	0.1 mL	407	Covance Core
MMS-570P-0100	Nestin Monoclonal Antibody	0.1 mL	298	Covance Core
MMS-570P-0200	Nestin Monoclonal Antibody	0.2 mL	498	Covance Core
MMS-580S-0200	GAPDH Monoclonal Antibody	0.2 mL	356	Covance Core
MMS-581S-0200	Fibrillarin Monoclonal Antibody	0.2 mL	356	Covance Core
MMS-582S-0200	Peripherin 8G2 Monoclonal Antibody	0.2 mL	356	Covance Core
MMS-583S-0200	Neurofilament NF-M Monoclonal Antibody	0.2 mL	356	Covance Core
MMS-584S-0200	Neurofilament NF-L Monoclonal Antibody	0.2 mL	356	Covance Core
MMS-585R-0100	Neurofilament NF-H Monoclonal Antibody	0.1 mL	356	Covance Core
MMS-586S-0200	Neurofilament α -Internexin/NF66 Monoclonal Antibody	0.2 mL	356	Covance Core
MMS-587S-0200	Neurofilament α -Internexin/NF66 Monoclonal Antibody	0.2 mL	356	Covance Core
MMS-588S-0200	GFAP Monoclonal Antibody	0.2 mL	356	Covance Core
MMS-589P-0200	Peripherin 7C5 Monoclonal Antibody	0.2 mL	356	Covance Core
MMS-604R-0100	Complement Receptor gC1q-R/p33 Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-606R-0100	Complement Receptor gC1q-R/p33 Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-610P-0250	Parathyroid Hormone Receptor, Opossum, Monoclonal Antibody, Purified	0.25 mL	467	Covance Core
MMS-612P-0500	Neuroendocrine (Chromogranin A) Monoclonal Antibody	0.5 mL	436	Covance Core
MMS-614R-0500	SNAP-25 (Synaptosomal-Associated Protein-25) Monoclonal Antibody	0.5 mL	486	Covance Core

CatalogNo	Product Name	Size	EUR	Product Line
MMS-616R-0500	Synaptobrevin Monoclonal Antibody	0.5 mL	486	Covance Core
MMS-618R-0500	Synaptophysin Monoclonal Antibody	0.5 mL	482	Covance Core
MMS-619R-0500	Syntaxin Monoclonal Antibody	0.5 mL	482	Covance Core
MMS-627R-0500	Neuronal Nicotinic Acetylcholine Receptor, α ;7-Subunit, Monoclonal Antibody	0.5 mL	467	Covance Core
MMS-645R-0100	ABCA3 (p180 Lamellar Body Protein) Monoclonal Antibody	0.1 mL	297	Covance Core
MMS-645R-0500	ABCA3 (p180 Lamellar Body Protein) Monoclonal Antibody	0.5 mL	552	Covance Core
MMS-660P-0250	Mucin (17B1) Monoclonal Antibody, Purified	0.25 mL	361	Covance Core
MMS-662P-0250	Mucin (17Q2) Monoclonal Antibody, Purified	0.25 mL	361	Covance Core
MPY-127R-0200	RNA Polymerase II Monoclonal Antibody Sampler	3 x 0.2 mL	752	Covance Core
MPY-661P-0500	Mucin, Matched Pair for ELISA	2 x 0.25 mL	600	Covance Core
MRB-435P-0100	Neuronal Class III β -Tubulin (TUJ1) Rabbit Monoclonal Antibody, Purified	0.1 mL	422	Covance Core
MRT-131P-0200	Maltose Binding Protein (MBP) Monoclonal Antibody, Purified	0.2 mL	400	Covance Core
MRT-204R-0200	PP2A Alpha Subunit (6F9) Monoclonal Antibody	0.2 mL	348	Covance Core
MRT-204R-0500	PP2A Alpha Subunit (6F9) Monoclonal Antibody	0.5 mL	482	Covance Core
MRT-207R-0200	PP2A Alpha Subunit (5H4) Monoclonal Antibody	0.2 mL	348	Covance Core
MRT-207R-0500	PP2A Alpha Subunit (5H4) Monoclonal Antibody	0.5 mL	482	Covance Core
MRT-208R-0200	PP2A Alpha Subunit (6G3) Monoclonal Antibody	0.2 mL	348	Covance Core
MRT-208R-0500	PP2A Alpha Subunit (6G3) Monoclonal Antibody	0.5 mL	482	Covance Core
MRT-601R-0500	Neuronal Nicotinic Acetylcholine Receptor, α ;8-Subunit, Monoclonal Antibody	0.5 mL	467	Covance Core
MRT-603R-0500	Neuronal Nicotinic Acetylcholine Receptor, α ;1, α ;3 and α ;5 Subunits,	0.5 mL	467	Covance Core
MRT-607R-0500	Neuronal Nicotinic Acetylcholine Receptor, α ;1-Subunit, Monoclonal Antibody	0.5 mL	467	Covance Core
MRT-609R-0500	Neuronal Nicotinic Acetylcholine Receptor, α ;1, α ;5 and α ;3 Subunits,	0.5 mL	467	Covance Core
MRT-611R-0500	Neuronal Nicotinic Acetylcholine Receptor α ;3-Subunits, Monoclonal Antibody	0.5 mL	467	Covance Core
MRT-613R-0500	Neuronal Nicotinic Acetylcholine Receptor α ;4-Subunits, Monoclonal Antibody	0.5 mL	467	Covance Core
MRT-623R-0500	Neuronal Nicotinic Acetylcholine Receptor α ;5-Subunits, Monoclonal Antibody	0.5 mL	467	Covance Core
MRT-629R-0500	Neuronal Nicotinic Acetylcholine Receptor, α ;7-Subunit, Monoclonal Antibody	0.5 mL	467	Covance Core
MRT-631R-0500	Neuronal Nicotinic Acetylcholine Receptor, α ;8-Subunits, Monoclonal Antibody	0.5 mL	467	Covance Core
MRT-633R-0500	Neuronal Nicotinic Acetylcholine Receptor β ;1-Subunits, Monoclonal Antibody	0.5 mL	467	Covance Core
MRT-639R-0500	Neuronal Nicotinic Acetylcholine Receptor, β ;2-Subunits, Monoclonal Antibody	0.5 mL	467	Covance Core
MRT-643R-0500	Neuronal Nicotinic Acetylcholine Receptor γ ;Subunits, Monoclonal Antibody	0.5 mL	467	Covance Core
PCK-153P-0100	Human Keratin 15 Polyclonal Antibody, Purified	0.1 mL	489	Covance Core
PCK-308P-0100	Neurofilament NF-L Polyclonal Antibody	0.1 mL	356	Covance Core
PCK-554P-0050	MAP2 Polyclonal Antibody, Purified	0.05 mL	422	Covance Core
PCK-591P-0100	GFAP Polyclonal Antibody, Purified	0.1 mL	549	Covance Core
PCK-592P-0050	Neurofilament NF-H Polyclonal Antibody, Purified	0.05 mL	422	Covance Core
PCK-593P-0100	Neurofilament NF-M Polyclonal Antibody, Purified	0.1 mL	356	Covance Core
PCK-594P-0100	Vimentin Polyclonal Antibody, Purified	0.1 mL	356	Covance Core
PEP-101P-1000	HA.11 Peptide, Purified	1.0 mg	251	Covance Core
PEP-101P-5	HA.11 Peptide, Purified	5.0 mg	482	Covance Core
PEP-115P-1000	GLU-GLU Peptide, Purified	1.0 mg	251	Covance Core
PEP-115P-5	GLU-GLU Peptide, Purified	5.0 mg	482	Covance Core
PEP-125P-1000	KT3 Peptide, Purified	1.0 mg	251	Covance Core
PEP-125P-5	KT3 Peptide, Purified	5.0 mg	482	Covance Core
PEP-130P-1000	AU1 Peptide, Purified	1.0 mg	251	Covance Core
PEP-130P-5	AU1 Peptide, Purified	5.0 mg	482	Covance Core
PEP-135P-1000	AU5 Peptide, Purified	1.0 mg	251	Covance Core
PEP-135P-5	AU5 Peptide, Purified	5.0 mg	482	Covance Core
PEP-150P-1000	9E10 (c-myc) Peptide, Purified	1.0 mg	251	Covance Core
PEP-150P-5	9E10 (c-myc) Peptide, Purified	5.0 mg	482	Covance Core
PEP-156P-1000	6-His Peptide, Purified	1.0 mg	251	Covance Core
PEP-156P-5	6-His Peptide, Purified	5.0 mg	482	Covance Core
PEP-435P-1000	Neuronal Class III β -Tubulin Peptide, Purified	1 mg	251	Covance Core
PGP-624C-0100	BACE1 C-terminal Guinea Pig Polyclonal Antibody	0.1 mL	467	Covance Core
PGT-180P-0200	cAMP (cyclic AMP), Polyclonal Antibody, Purified	0.2 mL	498	Covance Core
PGT-180P-0500	cAMP (cyclic AMP), Polyclonal Antibody, Purified	0.5 mL	752	Covance Core
PRB-101C-0200	HA.11 Polyclonal Antibody	0.2 mL	438	Covance Core
PRB-101C-0500	HA.11 Polyclonal Antibody	0.5 mL	660	Covance Core
PRB-101P-0200	HA.11 Polyclonal Antibody, Purified	0.2 mL	498	Covance Core
PRB-101P-0500	HA.11 Polyclonal Antibody, Purified	0.5 mL	752	Covance Core
PRB-105P-0200	Alpha B Crystallin Polyclonal Antibody, Purified	0.2 mL	467	Covance Core
PRB-105P-0500	Alpha B Crystallin Polyclonal Antibody, Purified	0.5 mL	630	Covance Core
PRB-106C-0200	Cre Recombinase Polyclonal Antibody	0.2 mL	438	Covance Core
PRB-112C-0200	GST Polyclonal Antibody	0.2 mL	438	Covance Core
PRB-113C-0100	Lamin A Polyclonal Antibody	0.1 mL	429	Covance Core
PRB-114C-0200	Giantin Polyclonal Antibody	0.2 mL	445	Covance Core
PRB-115C-0200	GLU-GLU Polyclonal Antibody	0.2 mL	438	Covance Core
PRB-117C-0100	Lamin C Polyclonal Antibody	0.1 mL	429	Covance Core
PRB-123C-0200	Eps15 (N-term), Polyclonal Antibody	0.2 mL	400	Covance Core
PRB-124C-0200	Eps15 (C-term), Polyclonal Antibody	0.2 mL	400	Covance Core
PRB-130P-0100	AU1 Polyclonal Antibody, Purified	0.1 mL	374	Covance Core
PRB-132C-0200	DYKDDDDK Polyclonal Antibody	0.2 mL	422	Covance Core
PRB-132P-0200	DYKDDDDK Polyclonal Antibody, Purified	0.2 mL	480	Covance Core
PRB-135C-0200	AU5 Polyclonal Antibody	0.2 mL	438	Covance Core
PRB-140C-0200	Involucrin Polyclonal Antibody	0.2 mL	422	Covance Core
PRB-144C-0200	GPP130 Polyclonal Antibody	0.2 mL	445	Covance Core
PRB-145P-0100	Loricrin Polyclonal Antibody, Purified	0.1 mL	489	Covance Core
PRB-149P-0100	Human Keratin 1 Polyclonal Antibody, Purified	0.1 mL	489	Covance Core

CatalogNo	Product Name	Size	EUR	Product Line
PRB-150C-0200	c-myc Polyclonal Antibody	0.2 mL	438	Covance Core
PRB-150C-0500	c-myc Polyclonal Antibody	0.5 mL	660	Covance Core
PRB-150P-0200	c-myc Polyclonal Antibody, Purified	0.2 mL	498	Covance Core
PRB-150P-0500	c-myc Polyclonal Antibody, Purified	0.5 mL	752	Covance Core
PRB-155P-0100	Mouse Keratin 14 (AF 64) Polyclonal Antibody, Purified	0.1 mL	489	Covance Core
PRB-156C-0200	6-His Polyclonal Antibody	0.2 mL	438	Covance Core
PRB-156C-0500	6-His Polyclonal Antibody	0.5 mL	660	Covance Core
PRB-159P-0100	Mouse Keratin 10 Polyclonal Antibody, Purified	0.1 mL	489	Covance Core
PRB-160P-0100	Mouse Keratin 5 (AF 138) Polyclonal Antibody, Purified	0.1 mL	489	Covance Core
PRB-163P-0100	Anti-Fluorescein Isothiocyanate (FITC) Polyclonal Antibody	0.1 mL	330	Covance Core
PRB-165P-0100	Mouse Keratin 1 (AF 109) Polyclonal Antibody, Purified	0.1 mL	489	Covance Core
PRB-167P-0100	Anti-Biotin Polyclonal Antibody	0.1 mL	330	Covance Core
PRB-169P-0100	Mouse Keratin 6 Polyclonal Antibody, Purified	0.1 mL	489	Covance Core
PRB-186P-0100	T7 Polyclonal Antibody	0.1 mL	374	Covance Core
PRB-189P-0100	V5 Polyclonal Antibody	0.1 mL	374	Covance Core
PRB-192P-0100	VSV-G Polyclonal Antibody	0.1 mL	374	Covance Core
PRB-212C-0200	Ran Polyclonal Antibody	0.2 mL	422	Covance Core
PRB-230C-0200	PIT-1 Polyclonal Antibody	0.2 mL	400	Covance Core
PRB-231P-0100	Hoxb-1 Polyclonal Antibody, Purified	0.1 mL	438	Covance Core
PRB-236P-0100	Krox-20 (Egr2) Polyclonal Antibody, Purified	0.1 mL	459	Covance Core
PRB-238C-0200	Prox1 Polyclonal Antibody	0.2 mL	452	Covance Core
PRB-250C-0200	KIN-28 Polyclonal Antibody	0.2 mL	400	Covance Core
PRB-252C-0200	Focal Adhesion Kinase (pp125FAK) Polyclonal Antibody	0.2 mL	400	Covance Core
PRB-255C-0200	GAL4 Polyclonal Antibody	0.2 mL	400	Covance Core
PRB-268C-0100	Ubiquitin Polyclonal Antibody	0.1 mL	356	Covance Core
PRB-271C-0100	UBE1L N-Terminal Polyclonal Antibody	0.1 mL	400	Covance Core
PRB-272C-0100	UBE1L C-Terminal Polyclonal Antibody	0.1 mL	400	Covance Core
PRB-276P-0200	Pax-2 Polyclonal Antibody, Purified	0.2 mL	498	Covance Core
PRB-278P-0100	Pax-6 Polyclonal Antibody, Purified	0.1 mL	459	Covance Core
PRB-282C-0100	Insulin Degrading Enzyme (IDE) Polyclonal Antibody	0.1 mL	400	Covance Core
PRB-291C-0200	Drosophila HP1 Polyclonal Antibody	0.2 mL	422	Covance Core
PRB-293C-0200	Drosophila HP1 Polyclonal Antibody	0.2 mL	422	Covance Core
PRB-297C-0200	Drosophila HP1 Polyclonal Antibody	0.2 mL	422	Covance Core
PRB-313C-0100	Human Nestin Repeat Polyclonal Antibody	0.1 mL	429	Covance Core
PRB-314C-0100	Mouse Nestin Repeat Polyclonal Antibody	0.1 mL	429	Covance Core
PRB-315C-0100	Mouse Nestin Tail Polyclonal Antibody	0.1 mL	429	Covance Core
PRB-317P-0100	Dopamine Decarboxylase Polyclonal Antibody	0.1 mL	429	Covance Core
PRB-319P-0100	Metabotropic Glutamate Receptor 1a Polyclonal Antibody	0.1 mL	429	Covance Core
PRB-320P-0100	Metabotropic Glutamate Receptor 2/3 Polyclonal Antibody	0.1 mL	429	Covance Core
PRB-321P-0100	Metabotropic Glutamate Receptor 5/1a Polyclonal Antibody	0.1 mL	429	Covance Core
PRB-323P-0100	Human DOPA Decarboxylase Polyclonal Antibody	0.1 mL	429	Covance Core
PRB-324P-0100	Dopamine Transporter, C-Terminus Polyclonal Antibody	0.1 mL	429	Covance Core
PRB-328P-0200	Glycine Receptor Polyclonal Antibody	0.2 mg	429	Covance Core
PRB-329P-0100	Vesicular Monoamine Transporter 2, C-Terminus Polyclonal Antibody	0.1 mL	429	Covance Core
PRB-330P-0100	Dopamine Transporter EL2 Polyclonal Antibody	0.1 mL	429	Covance Core
PRB-331P-0100	Voltage Gated Potassium Channel, Kv2.2 Subunit Polyclonal Antibody	0.1 mL	429	Covance Core
PRB-336C-0100	Human Nestin Tail Polyclonal Antibody	0.1 mL	429	Covance Core
PRB-354P-0100	Presenilin 1/N-Terminal (PS1/N-Term) Polyclonal Antibody, Purified	0.1 mL	422	Covance Core
PRB-355P-0100	Presenilin 1/C-Terminal (PS1/C-Term) Polyclonal Antibody, Purified	0.1 mL	422	Covance Core
PRB-362C-0100	ER81 Polyclonal Antibody	0.1 mL	429	Covance Core
PRB-364P-0100	Nicastrin N-Terminal (Nicastrin/N-Term) Polyclonal Antibody, Purified	0.1 mL	422	Covance Core
PRB-365P-0100	Nicastrin C-Terminal (Nicastrin/C-Term) Polyclonal Antibody, Purified	0.1 mL	422	Covance Core
PRB-417P-0100	Filaggrin Polyclonal Antibody, Purified	0.1 mL	489	Covance Core
PRB-432C-0200	Pericentrin Polyclonal Antibody	0.2 mL	422	Covance Core
PRB-433C-0200	γ-Tubulin Polyclonal Antibody	0.2 mL	422	Covance Core
PRB-435P-0100	Neuronal Class III β-Tubulin Polyclonal Antibody, Purified	0.1 mL	482	Covance Core
PRB-439P-0100	Nesprin-1 Polyclonal Antibody	0.1 mL	429	Covance Core
PRB-440P-0100	Nonmuscle Myosin II Heavy Chain A (MHC-A) Polyclonal Antibody, Purified	0.1 mL	467	Covance Core
PRB-445P-0100	Nonmuscle Myosin II Heavy Chain B (MHC-B) Polyclonal Antibody, Purified	0.1 mL	467	Covance Core
PRB-451P-0100	mAKAP Polyclonal Antibody	0.1 mL	429	Covance Core
PRB-452C-0200	MAD2 Polyclonal Antibody	0.2 mL	452	Covance Core
PRB-457C-0100	Myosin Phosphatase Polyclonal Antibody	0.1 mL	400	Covance Core
PRB-462P-050	Amyloid-β 1-40 Polyclonal Antibody, Purified	lyophilized	651	Covance Core
PRB-465P-050	Amyloid-β Polyclonal Antibody, Purified	lyophilized	651	Covance Core
PRB-497P-0250	S100A4 Polyclonal Antibody, Purified	0.25 mL	434	Covance Core
PRB-498C-0100	Amyloid-β 1-40 Polyclonal Antibody	0.1 mL	450	Covance Core
PRB-499C-0100	Amyloid-β 1-42 Polyclonal Antibody	0.1 mL	450	Covance Core
PRB-502P-0100	Tyrosine Hydroxylase [Ser40] Polyclonal Antibody, Phospho-Specific	0.1 mL	436	Covance Core
PRB-504P-0100	MEK 1/2 [Ser 218/222] Polyclonal Antibody, Phospho-Specific	0.1 mL	436	Covance Core
PRB-505P-0100	MARCKS [Ser152/156] Polyclonal Antibody, Phospho-specific	0.1 mL	436	Covance Core
PRB-506P-0100	Calcium/Calmodulin Kinase II [Thr286] Polyclonal Antibody, Phospho-Specific	0.1 mL	436	Covance Core
PRB-507P-0150	GluR1 [Ser831] Polyclonal Antibody, Phospho-specific	0.15 mL	436	Covance Core
PRB-508P-0100	ERK [p44/42 MAP Kinase] [Thr202/Tyr204] Polyclonal Antibody, Phospho-Specific	0.1 mL	436	Covance Core
PRB-509P-0150	GluR1 [Ser845] Polyclonal Antibody, Phospho-Specific	0.15 mL	436	Covance Core
PRB-510P-0100	p38 MAPK [Thr180/Tyr182] Polyclonal Antibody, Phospho-specific	0.1 mL	436	Covance Core
PRB-511P-0100	p53 [Ser392] Polyclonal Antibody, Phospho-specific	0.1 mL	436	Covance Core
PRB-512P-010	NMDA NR2B Polyclonal Antibody, Purified	0.01 mg	436	Covance Core

CatalogNo	Product Name	Size	EUR	Product Line
PRB-513P-010	NMDA NR2A Polyclonal Antibody, Purified	0.01 mg	436	Covance Core
PRB-514P-010	NMDA NR2C Polyclonal Antibody, Purified	0.01 mg	436	Covance Core
PRB-515P-0100	Tyrosine Hydroxylase Polyclonal Antibody	0.1 mL	436	Covance Core
PRB-516P-010	Synapsin I Polyclonal Antibody, Purified	0.01 mg	436	Covance Core
PRB-547C-0100	MAP2 Polyclonal Antibody	0.1 mL	422	Covance Core
PRB-548P-0100	GABA_A Receptor α4 Subunit N-terminus Polyclonal Antibody	0.1 mL	400	Covance Core
PRB-549P-0100	GABA_A Receptor α4 Subunit Cytosolic Loop Polyclonal Antibody	0.1 mL	400	Covance Core
PRB-550P-0100	Aph-1aL/C Terminal (245-265) Polyclonal Antibody, Purified	0.1 mL	498	Covance Core
PRB-551P-0100	Aph-1aL/S Loop (92-115) Polyclonal Antibody, Purified	0.1 mL	498	Covance Core
PRB-552P-0100	Aph-1b/C Terminal (245-257) Polyclonal Antibody, Purified	0.1 mL	498	Covance Core
PRB-553P-0100	Aph-1b/Loop Polyclonal Antibody, Purified	0.1 mL	498	Covance Core
PRB-558P-0100	GABA_A Receptor α6 Subunit C-Terminal Polyclonal Antibody	0.1 mL	467	Covance Core
PRB-559C-0050	GABA_A Receptor α6 Subunit Polyclonal Antibody	0.05 mL	400	Covance Core
PRB-560C-0100	PEN2 (N-term) Polyclonal Antibody	0.1 mL	422	Covance Core
PRB-562C-0050	GABA_A Receptor α2 Subunit Polyclonal Antibody	0.05 mL	400	Covance Core
PRB-563P-0100	GABA_A Receptor α1 Subunit N-Terminus Polyclonal Antibody	0.1 mL	467	Covance Core
PRB-564P-0100	GABA_A Receptor α1 Subunit C-Terminus Polyclonal Antibody	0.1 mL	467	Covance Core
PRB-567C-0050	GABA_A Receptor α1 Subunit Polyclonal Antibody	0.05 mL	400	Covance Core
PRB-568P-0100	GABA_A Receptor γ2 Subunit C-Terminus Loop Polyclonal Antit	0.1 mL	467	Covance Core
PRB-569C-0050	GABA_A Receptor γ2 Subunit Polyclonal Antibody	0.05 mL	400	Covance Core
PRB-570C-0100	Nestin Polyclonal Antibody	0.1 mL	429	Covance Core
PRB-571C-0100	GFAP Polyclonal Antibody	0.1 mL	356	Covance Core
PRB-572C-0100	Neurofilament α-Internexin/NF66 Polyclonal Antibody	0.1 mL	356	Covance Core
PRB-573C-0100	Neurofilament NF-H Polyclonal Antibody	0.1 mL	356	Covance Core
PRB-574C-0100	Neurofilament NF-L Polyclonal Antibody	0.1 mL	356	Covance Core
PRB-575C-0100	Neurofilament NF-M Polyclonal Antibody	0.1 mL	356	Covance Core
PRB-576C-0100	Panspecific Peripherin Polyclonal Antibody	0.1 mL	356	Covance Core
PRB-595P-0100	GABA_A Receptor β3 Subunit C-Terminus Polyclonal Antibody	0.1 mL	467	Covance Core
PRB-596P-0100	GABA_A Receptor β2 Subunit C-Terminus Polyclonal Antibody	0.1 mL	467	Covance Core
PRB-599C-0100	PEN2 (C-term) Polyclonal Antibody	0.1 mL	422	Covance Core
PRB-602C-0200	p75 Neurotrophin Receptor Polyclonal Antibody	0.2 mL	400	Covance Core
PRB-608P-0100	p75 Neurotrophin Receptor Polyclonal Antibody, Purified	0.1 mL	467	Covance Core
PRB-617C-0100	BACE1 C-terminal Rabbit Polyclonal Antibody	0.1 mL	467	Covance Core
PRB-620P-0100	Parathyroid Hormone Receptor Polyclonal Antibody, Peptide II	0.1 mL	467	Covance Core
PRB-622C-0100	ProBACE1 Polyclonal Antibody	0.1 mL	467	Covance Core
PRB-625P-0100	Parathyroid Hormone Receptor Polyclonal Antibody, Peptide I	0.1 mL	467	Covance Core
PRB-630P-0100	Parathyroid Hormone Receptor Polyclonal Antibody, Peptide IV	0.1 mL	467	Covance Core
PRB-635P-0100	Parathyroid Hormone (PTH) Receptor Polyclonal Antibody, Peptide VII	0.1 mL	467	Covance Core
PRB-640P-0100	Parathyroid Hormone Receptor, Human, Polyclonal Antibody	0.1 mL	467	Covance Core
PRO-133P-0050	Posi-Tag Epitope Tag Control Protein	0.05 mL	166	Covance Core
PSH-131P-0200	Maltose Binding Protein (MBP) Polyclonal Antibody, Purified	0.2 mL	498	Covance Core
PSH-131P-0500	Maltose Binding Protein (MBP) Polyclonal Antibody, Purified	0.5 mL	660	Covance Core
PSH-322P-0100	DOPA Decarboxylase Polyclonal Antibody	0.1 mL	429	Covance Core
PSH-325P-0100	Dopamine β-Hydroxylase, N-Terminus Polyclonal Antibody	0.1 mL	429	Covance Core
PSH-326P-0100	Dopamine β-Hydroxylase, C-Terminus Polyclonal Antibody	0.1 mL	429	Covance Core
PSH-327P-0100	Tryptophan Hydroxylase Polyclonal Antibody	0.1 mL	429	Covance Core
rPEP-579P-1000	Amyloid-β; Recombinant Peptide (1-40)	1.0 mg	400	Covance Core
rPEP-590P-1000	Amyloid-β; Recombinant Peptide (1-42)	1.0 mg	512	Covance Core
rPEP-597P-0500	Amyloid-β; Scrambled Recombinant Peptide (1-40)	0.5 mg	289	Covance Core
rPEP-598P-0500	Amyloid-β; Scrambled Recombinant Peptide (1-42)	0.5 mg	400	Covance Core
SIG-31010-5	PAMM (Permanent Aqueous Mounting Medium), 5 mL (SIGNET)	5.0 mL	55	IHC
SIG-31012-15	PAMM (Permanent Aqueous Mounting Medium), 15 mL (SIGNET)	15 mL	61	IHC
SIG-31015-5	MAYER'S Hematoxylin, 5 mL (SIGNET)	5.0 mL	54	IHC
SIG-31018-15	MAYER'S Hematoxylin, 15 mL (SIGNET)	15 mL	61	IHC
SIG-31020-1	PBS, 10x Concentrate (SIGNET)	1.0 L	161	IHC
SIG-31022-1000	AEC Chromogen Concentrate (SIGNET)	1.0 mL	105	IHC
SIG-31023-2000	AEC Chromogen Concentrate (SIGNET)	2.0 mL	101	IHC
SIG-31024-kit	AEC Chromogen System, 300 Tests (SIGNET)	kit	132	IHC
SIG-31026-kit	AEC Chromogen System, 1000 Tests (SIGNET)	kit	192	IHC
SIG-31027-3	AEC Chromogen Concentrate (SIGNET)	3.0 mL	132	IHC
SIG-31030-0100	AEC Substrate Buffer (SIGNET)	2 x 50 mL	82	IHC
SIG-31035-1	Staining Tray (Humidity Chamber) (SIGNET)	1 pk	61	IHC
SIG-31036-2	Staining Tray (Humidity Chamber)	2 pkg	87	IHC
SIG-31040-kit	DAB Chromogen System, 300 Test (SIGNET)	kit	141	IHC
SIG-31041-4	DAB Chromogen Concentrate (SIGNET)	4.0 mL	101	IHC
SIG-31042-kit	DAB Chromogen System, 1000 Test (SIGNET)	kit	192	IHC
SIG-31043-6	DAB Chromogen Concentrate (SIGNET)	6.0 mL	138	IHC
SIG-31044-15	DAB Substrate Buffer (SIGNET)	15 mL	83	IHC
SIG-31045-0100	DAB Substrate Buffer (SIGNET)	2 x 50 mL	86	IHC
SIG-31047-3	Boost DAB Enhancer, 50x (SIGNET)	3.0 mL	82	IHC
SIG-31048-kit	DAB 2 Component with Stabilizer (SIGNET)	kit	138	IHC
SIG-31049-0100	DAB Substrate Buffer with Stabilizer (SIGNET)	2 x 50 mL	109	IHC
SIG-31070-kit	FAST RED Chromogen System (SIGNET)	kit	152	IHC
SIG-31072-kit	FAST RED Chromogen System (SIGNET)	kit	192	IHC
SIG-31073-20	FAST RED Chromogen (SIGNET)	20 Tablets	152	IHC
SIG-31076-0100	FAST RED Substrate Buffer (SIGNET)	2 x 50 mL	86	IHC
SIG-31081-1	NBT/BCIP Chromogen System (SIGNET)	1 L	1346	IHC

CatalogNo	Product Name	Size	EUR	Product Line
SIG-31082-30	NBT/BCIP Chromogen System (SIGNET)	2 x 15 mL	152	IHC
SIG-31084-0100	NBT/BCIP Chromogen System (SIGNET)	2 x 50 mL	219	IHC
SIG-31095-kit	Xylene Resistant (X/R) Red System, 300 Tests (SIGNET)	kit	201	IHC
SIG-31097-kit	Xylene Resistant (X/R) Red System, 1000 Tests (SIGNET)	kit	522	IHC
SIG-3110-1000	α-1-Fetoprotein Polyclonal Antibody, Purified (SIGNET)	1.0 mL	234	IHC
SIG-3110-16	α-1-Fetoprotein Polyclonal Antibody, Purified (SIGNET)	L1 6mL	198	IHC
SIG-3110-26	α-1-Fetoprotein Polyclonal Antibody, Purified (SIGNET)	L2 6mL	176	IHC
SIG-3111-1000	Carcinoembryonic Antigen (CEA) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	438	IHC
SIG-3111-16	Carcinoembryonic Antigen (CEA) Monoclonal Antibody, Purified (SIGNET)	L1 6mL	234	IHC
SIG-3111-26	Carcinoembryonic Antigen (CEA) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	208	IHC
SIG-31120-50	Antibody Diluent Buffer (SIGNET)	50 mL	72	IHC
SIG-3112-1000	Carcinoembryonic Antigen (CEA) Polyclonal Antibody, Purified (SIGNET)	1.0 mL	234	IHC
SIG-31121-250	Antibody Diluent Buffer (SIGNET)	250 mL	215	IHC
SIG-3112-16	Carcinoembryonic Antigen (CEA) Polyclonal Antibody, Purified (SIGNET)	L1 6mL	193	IHC
SIG-3112-26	Carcinoembryonic Antigen (CEA) Polyclonal Antibody, Purified (SIGNET)	L2 6mL	173	IHC
SIG-3112-36	Carcinoembryonic Antigen (CEA) Polyclonal Antibody, Purified (SIGNET)	LV 6mL	284	IHC
SIG-31126-kit	Avidin/Biotin Blocking System (SIGNET)	kit	134	IHC
SIG-31128-15	Peroxide Blocking Reagent, 15mL (SIGNET)	15 mL	61	IHC
SIG-31130-50	Peroxide Blocking Reagent, 50mL (SIGNET)	50 mL	73	IHC
SIG-3113-1000	Epithelial Membrane Antigen (EMA) Monoclonal Antibody (SIGNET)	1.0 mL	289	IHC
SIG-3113-16	Epithelial Membrane Antigen (EMA) Monoclonal Antibody (SIGNET)	L1 6mL	272	IHC
SIG-3113-26	Epithelial Membrane Antigen (EMA) Monoclonal Antibody (SIGNET)	L2 6mL	234	IHC
SIG-3113-36	Epithelial Membrane Antigen (EMA) Monoclonal Antibody (SIGNET)	LV 6mL	284	IHC
SIG-3114-1000	Factor VIII Related Antigen Monoclonal Antibody (SIGNET)	1.0 mL	272	IHC
SIG-3114-16	Factor VIII Related Antigen Monoclonal Antibody (SIGNET)	L1 6mL	235	IHC
SIG-3114-26	Factor VIII Related Antigen Monoclonal Antibody (SIGNET)	L2 6mL	217	IHC
SIG-3115-1000	Factor VIII Related Antigen Polyclonal Antibody, Purified (SIGNET)	1.0 mL	242	IHC
SIG-3115-16	Factor VIII Related Antigen Polyclonal Antibody, Purified (SIGNET)	L1 6mL	217	IHC
SIG-3115-26	Factor VIII Related Antigen Polyclonal Antibody, Purified (SIGNET)	L2 6mL	201	IHC
SIG-3116-1000	HMB-45 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	572	IHC
SIG-3116-16	HMB-45 Monoclonal Antibody, Purified (SIGNET)	L1 6mL	335	IHC
SIG-3116-26	HMB-45 Monoclonal Antibody, Purified (SIGNET)	L2 6mL	301	IHC
SIG-31170-6	Normal Serum Block, 6 mL (SIGNET)	6.0 mL	66	IHC
SIG-3117-1000	Prostate Specific Antigen (PSA) Polyclonal Antibody, Purified (SIGNET)	1.0 mL	234	IHC
SIG-31171-15	Normal Serum Block, 15 mL (SIGNET)	15 mL	88	IHC
SIG-3117-16	Prostate Specific Antigen (PSA) Polyclonal Antibody, Purified (SIGNET)	L1 6mL	187	IHC
SIG-31172-50	Normal Serum Block, 50 mL (SIGNET)	50 mL	146	IHC
SIG-3117-26	Prostate Specific Antigen (PSA) Polyclonal Antibody, Purified (SIGNET)	L2 6mL	169	IHC
SIG-3117-36	Prostate Specific Antigen (PSA) Polyclonal Antibody, Purified (SIGNET)	LV 6mL	272	IHC
SIG-3118-26	S100 Monoclonal Antibody, Purified (SIGNET)	L2 6mL	294	IHC
SIG-3118-36	S100 Monoclonal Antibody, Purified (SIGNET)	LV 6mL	272	IHC
SIG-3119-1000	S100 Polyclonal Antibody, Purified (SIGNET)	1.0 mL	283	IHC
SIG-3119-16	S100 Polyclonal Antibody, Purified (SIGNET)	L1 6mL	215	IHC
SIG-3119-26	S100 Polyclonal Antibody, Purified (SIGNET)	L2 6mL	201	IHC
SIG-3120-1000	Myoglobin Polyclonal Antibody, Purified (SIGNET)	1.0 mL	223	IHC
SIG-3120-16	Myoglobin Polyclonal Antibody, Purified (SIGNET)	L1 6mL	206	IHC
SIG-3120-26	Myoglobin Polyclonal Antibody, Purified (SIGNET)	L2 6mL	207	IHC
SIG-3121-1000	Prostate Specific Antigen (PSA) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	283	IHC
SIG-3121-16	Prostate Specific Antigen (PSA) Monoclonal Antibody, Purified (SIGNET)	L1 6mL	234	IHC
SIG-3121-26	Prostate Specific Antigen (PSA) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	106	IHC
SIG-31900-50	Retrieve-All Antigen Unmasking System 3: Acidic, 100x (SIGNET)	50 mL	398	IHC
SIG-31902-0500	Retrieve-All Antigen Unmasking System 3: Acidic, 1x (SIGNET)	500 mL	91	IHC
SIG-31904-0500	Retrieve-All Antigen Unmasking System 3: Acidic, 10x (SIGNET)	500 mL	398	IHC
SIG-31910-50	Retrieve-All Antigen Unmasking System 1: Universal, 100x (SIGNET)	50 mL	383	IHC
SIG-31912-0500	Retrieve-All Antigen Unmasking System 1: Universal, 1x (SIGNET)	500 mL	74	IHC
SIG-31914-0500	Retrieve-All Antigen Unmasking System 1: Universal, 10x (SIGNET)	500 mL	383	IHC
SIG-31920-50	Retrieve-All Antigen Unmasking System 2: Basic, 100x (SIGNET)	50 mL	398	IHC
SIG-31922-0500	Retrieve-All Antigen Unmasking System 2: Basic, 1x (SIGNET)	500 mL	91	IHC
SIG-31924-0500	Retrieve-All Antigen Unmasking System 2: Basic, 10x (SIGNET)	500 mL	398	IHC
SIG-31932-0500	Sodium Citrate H.I.E.R., 1x (SIGNET)	500 mL	77	IHC
SIG-31934-0500	Sodium Citrate H.I.E.R., 10x (SIGNET)	500 mL	398	IHC
SIG-3210-1000	Adrenocorticotrophic Hormone Polyclonal Antibody (SIGNET)	1.0 mL	306	IHC
SIG-3210-16	Adrenocorticotrophic Hormone Polyclonal Antibody (SIGNET)	L1 6mL	224	IHC
SIG-3210-26	Adrenocorticotrophic Hormone Polyclonal Antibody (SIGNET)	L2 6mL	207	IHC
SIG-3211-1000	Calcitonin Polyclonal Antibody, Purified (SIGNET)	1.0 mL	223	IHC
SIG-3211-16	Calcitonin Polyclonal Antibody, Purified (SIGNET)	L1 6mL	221	IHC
SIG-3211-26	Calcitonin Polyclonal Antibody, Purified (SIGNET)	L2 6mL	185	IHC
SIG-3213-1000	Human Chorionic Gonadotropin-β Polyclonal Antibody, Purified (SIGNET)	1.0 mL	223	IHC
SIG-3213-16	Human Chorionic Gonadotropin-β Polyclonal Antibody, Purified (SIGNET)	L1 6mL	221	IHC
SIG-3213-26	Human Chorionic Gonadotropin-β Polyclonal Antibody, Purified (SIGNET)	L2 6mL	185	IHC
SIG-3214-1000	Human Chorionic Gonadotropin (hCG) Polyclonal Antibody, Purified (SIGNET)	1.0 mL	221	IHC
SIG-3214-16	Human Chorionic Gonadotropin (hCG) Polyclonal Antibody, Purified (SIGNET)	L1 6mL	174	IHC
SIG-3214-26	Human Chorionic Gonadotropin (hCG) Polyclonal Antibody, Purified (SIGNET)	L2 6mL	156	IHC
SIG-3215-1000	Chromogranin A Monoclonal Antibody, Purified (SIGNET)	1.0 mL	299	IHC
SIG-3215-16	Chromogranin A Monoclonal Antibody, Purified (SIGNET)	L1 6mL	283	IHC
SIG-3215-26	Chromogranin A Monoclonal Antibody, Purified (SIGNET)	L2 6mL	249	IHC
SIG-3215-36	Chromogranin A Monoclonal Antibody, Purified (SIGNET)	LV 6mL	284	IHC

CatalogNo	Product Name	Size	EUR	Product Line
SIG-3217-1000	FSH Polyclonal Antibody, Purified (SIGNET)	1.0 mL	290	IHC
SIG-3217-16	FSH Polyclonal Antibody, Purified (SIGNET)	L1 6mL	228	IHC
SIG-3217-26	FSH Polyclonal Antibody, Purified (SIGNET)	L2 6mL	223	IHC
SIG-3219-1000	Glucagon Polyclonal Antibody, Purified (SIGNET)	1.0 mL	290	IHC
SIG-3219-16	Glucagon Polyclonal Antibody, Purified (SIGNET)	L1 6mL	228	IHC
SIG-3219-26	Glucagon Polyclonal Antibody, Purified (SIGNET)	L2 6mL	223	IHC
SIG-3220-1000	Growth Hormone (GH) Polyclonal Antibody, Purified (SIGNET)	1.0 mL	290	IHC
SIG-3220-16	Growth Hormone (GH) Polyclonal Antibody, Purified (SIGNET)	L1 6mL	228	IHC
SIG-3220-26	Growth Hormone (GH) Polyclonal Antibody, Purified (SIGNET)	L2 6mL	223	IHC
SIG-32230-kit	Ultrasreptavidin (USA) Horseradish Peroxidase (HRP) 50 Test, Rabbit, AEC Detection kit		258	IHC
SIG-32232-kit	Ultrasreptavidin (USA) Horseradish Peroxidase (HRP) 50 Test, Rabbit, DAB Detection kit		258	IHC
SIG-32234-kit	Ultrasreptavidin (USA) Horseradish Peroxidase (HRP) 150 Test, Rabbit, Detection Kit kit		398	IHC
SIG-32236-kit	Ultrasreptavidin (USA) Horseradish Peroxidase (HRP) 500 Test, Rabbit, Detection Kit kit		613	IHC
SIG-32238-kit	Ultrasreptavidin (USA) Horseradish Peroxidase (HRP) 50 Test, Murine, AEC Detection kit		258	IHC
SIG-32240-kit	Ultrasreptavidin (USA) Horseradish peroxidase (HRP) 50 Test, Murine, DAB Detection kit		258	IHC
SIG-3224-1000	Parathyroid Hormone (PTH) Polyclonal Antibody, Purified (SIGNET)	1.0 mL	316	IHC
SIG-3224-16	Parathyroid Hormone (PTH) Polyclonal Antibody, Purified (SIGNET)	L1 6mL	242	IHC
SIG-3224-26	Parathyroid Hormone (PTH) Polyclonal Antibody, Purified (SIGNET)	L2 6mL	217	IHC
SIG-32242-kit	Ultrasreptavidin (USA) Horseradish peroxidase (HRP) 150 Test, Murine, Detection Kit kit		398	IHC
SIG-32246-kit	Ultrasreptavidin (USA) Horseradish peroxidase (HRP) 500 Test, Murine, Detection Kit kit		613	IHC
SIG-32248-kit	Ultrasreptavidin (USA) Horseradish peroxidase (HRP) 50 Test, Multi-Species, AEC De kit		258	IHC
SIG-32250-kit	Ultrasreptavidin (USA) Horseradish peroxidase (HRP) 50 Test, Multi-Species, DAB De kit		258	IHC
SIG-3225-1000	Prolactin Polyclonal Antibody, Purified (SIGNET)	1.0 mL	316	IHC
SIG-3225-16	Prolactin Polyclonal Antibody, Purified (SIGNET)	L1 6mL	242	IHC
SIG-3225-26	Prolactin Polyclonal Antibody, Purified (SIGNET)	L2 6mL	217	IHC
SIG-32252-kit	Ultrasreptavidin (USA) Horseradish peroxidase (HRP) 150 Test, Multi-Species, Detecti kit		398	IHC
SIG-32254-kit	Ultrasreptavidin (USA) Horseradish peroxidase (HRP) 500 Test, Multi-Species, Detecti kit		589	IHC
SIG-32255-kit	Ultrasreptavidin (USA) Horseradish peroxidase (HRP) 1000 Test, Multi-Species, AEC I kit		1139	IHC
SIG-32256-kit	Ultrasreptavidin (USA) Horseradish peroxidase (HRP) 1000 Test, Multi-Species, DAB I kit		1139	IHC
SIG-3226-1000	Serotonin Polyclonal Antibody (SIGNET)	1.0 mL	316	IHC
SIG-3226-16	Serotonin Polyclonal Antibody (SIGNET)	L1 6mL	242	IHC
SIG-3226-26	Serotonin Polyclonal Antibody (SIGNET)	L2 6mL	217	IHC
SIG-3227-1000	Somatostatin Polyclonal Antibody (SIGNET)	1.0 mL	316	IHC
SIG-3227-16	Somatostatin Polyclonal Antibody (SIGNET)	L1 6mL	242	IHC
SIG-3227-26	Somatostatin Polyclonal Antibody (SIGNET)	L2 6mL	217	IHC
SIG-3228-1000	Thyroglobulin Polyclonal Antibody, Purified (SIGNET)	1.0 mL	316	IHC
SIG-3228-16	Thyroglobulin Polyclonal Antibody, Purified (SIGNET)	L1 6mL	242	IHC
SIG-3228-26	Thyroglobulin Polyclonal Antibody, Purified (SIGNET)	L2 6mL	217	IHC
SIG-3230-1000	Thyroid Stimulating Hormone (TSH) Polyclonal Antibody, Purified (SIGNET)	1.0 mL	316	IHC
SIG-3230-16	Thyroid Stimulating Hormone (TSH) Polyclonal Antibody, Purified (SIGNET)	L1 6mL	242	IHC
SIG-3230-26	Thyroid Stimulating Hormone (TSH) Polyclonal Antibody, Purified (SIGNET)	L2 6mL	217	IHC
SIG-32320-1	Capillary Action Buffer 1, 1L (SIGNET)	1.0 L	168	IHC
SIG-32330-1	Capillary Action Buffer 2/3, 1L (SIGNET)	1.0 L	168	IHC
SIG-32340-1	Capillary Action Water Wash, 1L (SIGNET)	1.0 L	168	IHC
SIG-32350-pkg	Capillary Action Trays (SIGNET)	250/pkg	482	IHC
SIG-32360-50	Capillary Action Absorbant Pads; 50 pcs (SIGNET)	50 pcs	412	New price April 2009
SIG-32380-0500	Tris-Buffered Saline 20X (1.0 M) w/ Tween 20 (SIGNET)	500 mL	148	IHC
SIG-32391-0500	Tris-Buffered Saline 20X (1.0 M) (SIGNET)	500 mL	150	IHC
SIG-32400-kit	Capillary Action Detection 700 Test HRP, AEC (SIGNET)	kit	1757	IHC
SIG-32402-kit	Capillary Action Detection 700 Test HRP, DAB (SIGNET)	700 test kit	1757	IHC
SIG-3250-1000	α-1-Antichymotrypsin Polyclonal Antibody, Purified (SIGNET)	1.0 mL	284	IHC
SIG-3250-16	α-1-Antichymotrypsin Polyclonal Antibody, Purified (SIGNET)	L1 6mL	221	IHC
SIG-3250-26	α-1-Antichymotrypsin Polyclonal Antibody, Purified (SIGNET)	L2 6mL	201	IHC
SIG-3251-1000	α-1-Antitrypsin Polyclonal Antibody, Purified (SIGNET)	1.0 mL	284	IHC
SIG-3251-16	α-1-Antitrypsin Polyclonal Antibody, Purified (SIGNET)	L1 6mL	221	IHC
SIG-3251-26	α-1-Antitrypsin Polyclonal Antibody, Purified (SIGNET)	L2 6mL	201	IHC
SIG-3255-1000	Lysozyme Polyclonal Antibody, Purified (SIGNET)	1.0 mL	284	IHC
SIG-3255-16	Lysozyme Polyclonal Antibody, Purified (SIGNET)	L1 6mL	221	IHC
SIG-3255-26	Lysozyme Polyclonal Antibody, Purified (SIGNET)	L2 6mL	201	IHC
SIG-3256-1000	Neuron Specific Enolase (NSE) Monoclonal Antibody (SIGNET)	1.0 mL	654	IHC
SIG-3256-16	Neuron Specific Enolase (NSE) Monoclonal Antibody (SIGNET)	L1 6mL	287	IHC
SIG-3256-26	Neuron Specific Enolase (NSE) Monoclonal Antibody (SIGNET)	L2 6mL	251	IHC
SIG-3257-1000	Neuron Specific Enolase (NSE) Polyclonal Antibody, Purified (SIGNET)	1.0 mL	439	IHC
SIG-3257-16	Neuron Specific Enolase (NSE) Polyclonal Antibody, Purified (SIGNET)	L1 6mL	379	IHC
SIG-3257-26	Neuron Specific Enolase (NSE) Polyclonal Antibody, Purified (SIGNET)	L2 6mL	280	IHC
SIG-3258-1000	Placental Alkaline Phosphatase (PLAP) Polyclonal Antibody, Purified (SIGNET)	1.0 mL	439	IHC
SIG-3258-16	Placental Alkaline Phosphatase (PLAP) Polyclonal Antibody, Purified (SIGNET)	L1 6mL	398	IHC
SIG-3258-26	Placental Alkaline Phosphatase (PLAP) Polyclonal Antibody, Purified (SIGNET)	L2 6mL	339	IHC
SIG-3259-1000	Prostatic Acid Phosphatase (PSAP) Polyclonal Antibody, Purified (SIGNET)	1.0 mL	243	IHC
SIG-3259-16	Prostatic Acid Phosphatase (PSAP) Polyclonal Antibody, Purified (SIGNET)	L1 6mL	217	IHC
SIG-3259-26	Prostatic Acid Phosphatase (PSAP) Polyclonal Antibody, Purified (SIGNET)	L2 6mL	192	IHC
SIG-32600-kit	ACUITY Biotin Free Polymer Detection HRP System 50 Test AEC (SIGNET)	kit	358	IHC
SIG-3260-26	Synaptophysin (SYP) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	438	IHC
SIG-32602-kit	ACUITY Biotin Free Polymer Detection HRP System 50 Test DAB (SIGNET)	kit	358	IHC
SIG-32604-kit	ACUITY Biotin Free Polymer Detection HRP System 150 Tests (SIGNET)	kit	687	IHC
SIG-32606-kit	ACUITY Biotin Free Polymer Detection HRP 500 Tests (SIGNET)	kit	1016	IHC
SIG-32610-kit	ACUITY Biotin Free Polymer Detection HRP (Cap Gap) 700 Tests (SIGNET)	kit	2168	IHC

CatalogNo	Product Name	Size	EUR	Product Line
SIG-3261-1000	Synaptophysin (SYP) Polyclonal Antibody, Purified (SIGNET)	1.0 mL	521	IHC
SIG-3261-26	Synaptophysin (SYP) Polyclonal Antibody, Purified (SIGNET)	L2 6mL	356	IHC
SIG-32612-kit	ACUITY Biotin Free Polymer Detection HRP (Cap Gap) Complete 700 Tests (SIGNET) kit		3323	IHC
SIG-32614-kit	ACUITY Biotin Free Polymer Detection HRP (Cap Gap) Complete 5,000 Tests (SIGNET) kit		25743	IHC
SIG-32702-kit	ACUITY Biotin Free Mouse-on-Mouse Polymer Detection System HRP 50 Test/DAB Kit kit		385	IHC
SIG-32704-kit	ACUITY Biotin Free Mouse-on-Mouse Polymer Detection System HRP 150 Test Kit	kit	978	IHC
SIG-32706-kit	ACUITY Biotin Free Mouse-on-Mouse Polymer Detection System HRP 500 Test Kit	kit	2799	IHC
SIG-32820-kit	ACUITYAdvanced Mouse-on-Mouse Polymer Detection System 50 Test DAB kit	kit	316	IHC
SIG-32822-kit	ACUITYAdvanced Mouse-on-Mouse Polymer Detection System 150 Test DAB kit	kit	796	IHC
SIG-32824-kit	ACUITYAdvanced Mouse-on-Mouse Polymer Detection System 500 Test DAB kit	kit	2268	IHC
SIG-32900-kit	ACUITYAdvanced Biotin Free Polymer Detection System 50 Test AEC kit	kit	294	IHC
SIG-32902-kit	ACUITYAdvanced Biotin Free Polymer Detection System 50 Test DAB kit	kit	294	IHC
SIG-32904-kit	ACUITYAdvanced Biotin Free Polymer Detection System 150 Test	kit	560	IHC
SIG-32906-kit	ACUITYAdvanced Biotin Free Polymer Detection System 500 Test	kit	827	IHC
SIG-32910-kit	ACUITYAdvanced Biotin Free Polymer Detection System (CAP GAP) 700 Test kit	kit	1759	IHC
SIG-32912-kit	ACUITYAdvanced Biotin Free Polymer Detection System (CAP GAP) Complete 700 Test kit	kit	2690	IHC
SIG-32914-kit	ACUITYAdvanced Biotin Free Polymer Detection System Complete 10,000 Test kit	kit	20810	IHC
SIG-3310-1000	BG-1 Monoclonal Antibody (SIGNET)	1.0 mL	448	IHC
SIG-3311-1000	BG-2 Monoclonal Antibody (SIGNET)	1.0 mL	448	IHC
SIG-3312-1000	BG-3 Monoclonal Antibody (SIGNET)	1.0 mL	448	IHC
SIG-3313-1000	BG-4 Monoclonal Antibody (SIGNET)	1.0 mL	448	IHC
SIG-3314-1000	BG-5 Monoclonal Antibody (SIGNET)	1.0 mL	448	IHC
SIG-3315-1000	BG-6 Monoclonal Antibody (SIGNET)	1.0 mL	448	IHC
SIG-3316-1000	BG-7 Monoclonal Antibody (SIGNET)	1.0 mL	448	IHC
SIG-3317-1000	BG-8 Monoclonal Antibody (SIGNET)	1.0 mL	448	IHC
SIG-3323-1000	CD15 Monoclonal Antibody (SIGNET)	1.0 mL	398	IHC
SIG-3323-16	CD15 Monoclonal Antibody (SIGNET)	L1 6mL	234	IHC
SIG-3323-26	CD15 Monoclonal Antibody (SIGNET)	L2 6mL	217	IHC
SIG-3324-1000	CD20 B Cell Monoclonal Antibody (SIGNET)	1.0 mL	398	IHC
SIG-3324-16	CD20 B Cell Monoclonal Antibody (SIGNET)	L1 6mL	234	IHC
SIG-3324-26	CD20 B Cell Monoclonal Antibody (SIGNET)	L2 6mL	217	IHC
SIG-3324-36	CD20 B Cell Monoclonal Antibody (SIGNET)	LV 6mL	275	IHC
SIG-3325-1000	CD30 Ki Antigen Monoclonal Antibody (SIGNET)	1.0 mL	582	IHC
SIG-3325-26	CD30 Ki Antigen Monoclonal Antibody (SIGNET)	L2 6mL	233	IHC
SIG-3326-1000	CD34 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	518	IHC
SIG-3326-26	CD34 Monoclonal Antibody, Purified (SIGNET)	L2 6mL	417	IHC
SIG-3327-1000	CD43 T-Cell Monoclonal Antibody (SIGNET)	1.0 mL	522	IHC
SIG-3327-16	CD43 T-Cell Monoclonal Antibody (SIGNET)	L1 6mL	439	IHC
SIG-3327-26	CD43 T-Cell Monoclonal Antibody (SIGNET)	L2 6mL	398	IHC
SIG-3327-36	CD43 T-Cell Monoclonal Antibody (SIGNET)	LV 6mL	439	IHC
SIG-3330-1000	Leukocyte Common Antigen Monoclonal Antibody (SIGNET)	1.0 mL	316	IHC
SIG-3330-16	Leukocyte Common Antigen Monoclonal Antibody (SIGNET)	L1 6mL	272	IHC
SIG-3330-26	Leukocyte Common Antigen Monoclonal Antibody (SIGNET)	L2 6mL	264	IHC
SIG-3330-36	Leukocyte Common Antigen Monoclonal Antibody (SIGNET)	LV 6mL	272	IHC
SIG-3331-1000	CD45RO Monoclonal Antibody (SIGNET)	1.0 mL	438	IHC
SIG-3331-16	CD45RO Monoclonal Antibody (SIGNET)	L1 6mL	352	IHC
SIG-3331-26	CD45RO Monoclonal Antibody (SIGNET)	L2 6mL	280	IHC
SIG-3331-36	CD45RO Monoclonal Antibody (SIGNET)	LV 6mL	238	IHC
SIG-3333-26	CDw75 (B Cell) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	275	IHC
SIG-3334-26	HLA-DR Monoclonal Antibody, Purified (SIGNET)	L2 6mL	275	IHC
SIG-3335-26	Macrophage Antigen Monoclonal Antibody (SIGNET)	L2 6mL	275	IHC
SIG-3352-26	Bcl-2 Oncoprotein Monoclonal Antibody (SIGNET)	L2 6mL	439	IHC
SIG-3410-1000	Immunoglobulin A (IgA) Polyclonal Antibody, Purified (SIGNET)	1.0 mL	188	IHC
SIG-3410-16	Immunoglobulin A (IgA) Polyclonal Antibody, Purified (SIGNET)	L1 6mL	153	IHC
SIG-3410-26	Immunoglobulin A (IgA) Polyclonal Antibody, Purified (SIGNET)	L2 6mL	132	IHC
SIG-3411-1000	Immunoglobulin G (IgG) Polyclonal Antibody, Purified (SIGNET)	1.0 mL	188	IHC
SIG-3411-16	Immunoglobulin G (IgG) Polyclonal Antibody, Purified (SIGNET)	L1 6mL	153	IHC
SIG-3411-26	Immunoglobulin G (IgG) Polyclonal Antibody, Purified (SIGNET)	L2 6mL	132	IHC
SIG-3414-1000	Immunoglobulin M (IgM) Polyclonal Antibody, Purified (SIGNET)	1.0 mL	188	IHC
SIG-3414-16	Immunoglobulin M (IgM) Polyclonal Antibody, Purified (SIGNET)	L1 6mL	153	IHC
SIG-3414-26	Immunoglobulin M (IgM) Polyclonal Antibody, Purified (SIGNET)	L2 6mL	132	IHC
SIG-3415-1000	Kappa Monoclonal Antibody (SIGNET)	1.0 mL	271	IHC
SIG-3415-16	Kappa Monoclonal Antibody (SIGNET)	L1 6mL	234	IHC
SIG-3415-26	Kappa Monoclonal Antibody (SIGNET)	L2 6mL	217	IHC
SIG-3416-1000	Kappa Polyclonal Antibody, Purified (SIGNET)	1.0 mL	175	IHC
SIG-3416-16	Kappa Polyclonal Antibody, Purified (SIGNET)	L1 6mL	152	IHC
SIG-3416-26	Kappa Polyclonal Antibody, Purified (SIGNET)	L2 6mL	142	IHC
SIG-3416-36	Kappa Polyclonal Antibody, Purified (SIGNET)	LV 6mL	275	IHC
SIG-3417-1000	Lambda Monoclonal Antibody (SIGNET)	1.0 mL	270	IHC
SIG-3417-16	Lambda Monoclonal Antibody (SIGNET)	L1 6mL	234	IHC
SIG-3417-26	Lambda Monoclonal Antibody (SIGNET)	L2 6mL	217	IHC
SIG-3418-1000	Lambda Polyclonal Antibody, Purified (SIGNET)	1.0 mL	175	IHC
SIG-3418-16	Lambda Polyclonal Antibody, Purified (SIGNET)	L1 6mL	152	IHC
SIG-3418-26	Lambda Polyclonal Antibody, Purified (SIGNET)	L2 6mL	142	IHC
SIG-3418-36	Lambda Polyclonal Antibody, Purified (SIGNET)	LV 6mL	275	IHC
SIG-3430-1000	Cytomegalovirus (CMV) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	472	IHC
SIG-3430-16	Cytomegalovirus (CMV) Monoclonal Antibody, Purified (SIGNET)	L1 6mL	356	IHC

CatalogNo	Product Name	Size	EUR	Product Line
SIG-3430-26	Cytomegalovirus (CMV) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	272	IHC
SIG-3431-26	H. pylori Polyclonal Antibody, Purified (SIGNET)	L2 6mL	272	IHC
SIG-3432-1000	Hepatitis B Core Antigen (HBcAg) Polyclonal Antibody, Purified (SIGNET)	1.0 mL	356	IHC
SIG-3432-16	Hepatitis B Core Antigen (HBcAg) Polyclonal Antibody, Purified (SIGNET)	L1 6mL	312	IHC
SIG-3432-26	Hepatitis B Core Antigen (HBcAg) Polyclonal Antibody, Purified (SIGNET)	L2 6mL	275	IHC
SIG-3433-1000	Hepatitis B Surface Antigen (HBsAg) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	366	IHC
SIG-3433-16	Hepatitis B Surface Antigen (HBsAg) Monoclonal Antibody, Purified (SIGNET)	L1 6mL	275	IHC
SIG-3433-26	Hepatitis B Surface Antigen (HBsAg) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	249	IHC
SIG-3434-1000	Hepatitis C Virus (HCV) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	921	IHC
SIG-3434-26	Hepatitis C Virus (HCV) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	518	IHC
SIG-3435-1000	Herpes Simplex Virus (HSV) Polyclonal Antibody, Purified (SIGNET)	1.0 mL	234	IHC
SIG-3435-16	Herpes Simplex Virus (HSV) Polyclonal Antibody, Purified (SIGNET)	L1 6mL	201	IHC
SIG-3435-26	Herpes Simplex Virus (HSV) Polyclonal Antibody, Purified (SIGNET)	L2 6mL	192	IHC
SIG-3436	N'est plus disponible			
SIG-3460-1000	α-Smooth Muscle Actin Monoclonal Antibody (SIGNET)	1.0 mL	358	IHC
SIG-3460-16	α-Smooth Muscle Actin Monoclonal Antibody (SIGNET)	L1 6mL	316	IHC
SIG-3460-26	α-Smooth Muscle Actin Monoclonal Antibody (SIGNET)	L2 6mL	292	IHC
SIG-3461-1000	Muscle Specific Actin Monoclonal Antibody (SIGNET)	1.0 mL	431	IHC
SIG-3461-16	Muscle Specific Actin Monoclonal Antibody (SIGNET)	L1 6mL	358	IHC
SIG-3461-26	Muscle Specific Actin Monoclonal Antibody (SIGNET)	L2 6mL	316	IHC
SIG-3462-1000	Cytokeratin (AE-1) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	275	IHC
SIG-3462-16	Cytokeratin (AE-1) Monoclonal Antibody, Purified (SIGNET)	L1 6mL	193	IHC
SIG-3462-26	Cytokeratin (AE-1) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	152	IHC
SIG-3463-1000	Cytokeratin (AE-3) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	275	IHC
SIG-3463-16	Cytokeratin (AE-3) Monoclonal Antibody, Purified (SIGNET)	L1 6mL	193	IHC
SIG-3463-26	Cytokeratin (AE-3) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	152	IHC
SIG-3464-1000	Pan-Cytokeratin (AE1/AE3) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	275	IHC
SIG-3464-16	Pan-Cytokeratin (AE1/AE3) Monoclonal Antibody, Purified (SIGNET)	L1 6mL	193	IHC
SIG-3464-26	Pan-Cytokeratin (AE1/AE3) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	152	IHC
SIG-3464-36	Pan-Cytokeratin (AE1/AE3) Monoclonal Antibody, Purified (SIGNET)	LV 6mL	275	IHC
SIG-3465-1000	Cytokeratin (OSCAR) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	482	IHC
SIG-3465-16	Cytokeratin (OSCAR) Monoclonal Antibody, Purified (SIGNET)	L1 6mL	325	IHC
SIG-3465-26	Cytokeratin (OSCAR) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	275	IHC
SIG-3469-1000	Desmin Monoclonal Antibody (SIGNET)	1.0 mL	383	IHC
SIG-3469-16	Desmin Monoclonal Antibody (SIGNET)	L1 6mL	290	IHC
SIG-3469-26	Desmin Monoclonal Antibody (SIGNET)	L2 6mL	275	IHC
SIG-3469-36	Desmin Monoclonal Antibody (SIGNET)	LV 6mL	292	IHC
SIG-3470-1000	GFAP (6F2) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	407	IHC
SIG-3470-16	GFAP (6F2) Monoclonal Antibody, Purified (SIGNET)	L1 6mL	275	IHC
SIG-3470-26	GFAP (6F2) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	234	IHC
SIG-3471-1000	Glial Fibrillary Acid Protein (GFAP) Polyclonal Antibody, Purified (SIGNET)	1.0 mL	316	IHC
SIG-3471-16	Glial Fibrillary Acid Protein (GFAP) Polyclonal Antibody, Purified (SIGNET)	L1 6mL	193	IHC
SIG-3471-26	Glial Fibrillary Acid Protein (GFAP) Polyclonal Antibody, Purified (SIGNET)	L2 6mL	175	IHC
SIG-3471-36	Glial Fibrillary Acid Protein (GFAP) Polyclonal Antibody, Purified (SIGNET)	LV 6mL	275	IHC
SIG-3472-1000	Myosin (MY-32) Monoclonal Antibody (SIGNET)	1.0 mL	383	IHC
SIG-3472-16	Myosin (MY-32) Monoclonal Antibody (SIGNET)	L1 6mL	234	IHC
SIG-3472-26	Myosin (MY-32) Monoclonal Antibody (SIGNET)	L2 6mL	217	IHC
SIG-3473-1000	Neurofilament (2F11) Monoclonal Antibody (SIGNET)	1.0 mL	383	IHC
SIG-3473-16	Neurofilament (2F11) Monoclonal Antibody (SIGNET)	L1 6mL	234	IHC
SIG-3473-26	Neurofilament (2F11) Monoclonal Antibody (SIGNET)	L2 6mL	275	IHC
SIG-3474-1000	Vimentin (V9) Monoclonal Antibody (SIGNET)	1.0 mL	383	IHC
SIG-3474-16	Vimentin (V9) Monoclonal Antibody (SIGNET)	L1 6mL	234	IHC
SIG-3474-26	Vimentin (V9) Monoclonal Antibody (SIGNET)	L2 6mL	275	IHC
SIG-3474-36	Vimentin (V9) Monoclonal Antibody (SIGNET)	LV 6mL	292	IHC
SIG-3512-1000	c-< >erb</> B2 Monoclonal Antibody (SIGNET)	1.0 mL	662	IHC
SIG-3512-16	c-< >erb</> B2 Monoclonal Antibody (SIGNET)	L1 6mL	518	IHC
SIG-3512-26	c-< >erb</> B2 Monoclonal Antibody (SIGNET)	L2 6mL	473	IHC
SIG-3513-1000	Epidermal Growth Factor Receptor (F4) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	568	IHC
SIG-3513-16	Epidermal Growth Factor Receptor (F4) Monoclonal Antibody, Purified (SIGNET)	L1 6mL	431	IHC
SIG-3513-26	Epidermal Growth Factor Receptor (F4) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	366	IHC
SIG-3515-26	Estrogen Receptor (1D5) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	582	IHC
SIG-3517-1000	p53 (BP53-12) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	635	IHC
SIG-3517-16	p53 (BP53-12) Monoclonal Antibody, Purified (SIGNET)	L1 6mL	603	IHC
SIG-3517-26	p53 (BP53-12) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	546	IHC
SIG-3518-1000	p53 (DO1) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	635	IHC
SIG-3518-16	p53 (DO1) Monoclonal Antibody, Purified (SIGNET)	L1 6mL	603	IHC
SIG-3518-26	p53 (DO1) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	546	IHC
SIG-3519-1000	p53 (DO7) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	635	IHC
SIG-3519-16	p53 (DO7) Monoclonal Antibody, Purified (SIGNET)	L1 6mL	603	IHC
SIG-3519-26	p53 (DO7) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	546	IHC
SIG-3520-1000	p53 CM1 Polyclonal Antibody (SIGNET)	1.0 mL	747	IHC
SIG-3520-16	p53 CM1 Polyclonal Antibody (SIGNET)	L1 6mL	546	IHC
SIG-3520-26	p53 CM1 Polyclonal Antibody (SIGNET)	L2 6mL	496	IHC
SIG-3521-26	Progesterone Receptor (PR) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	603	IHC
SIG-3523-1000	Proliferating Cell Nuclear Antigen (PCNA) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	493	IHC
SIG-3523-16	Proliferating Cell Nuclear Antigen (PCNA) Monoclonal Antibody, Purified (SIGNET)	L1 6mL	316	IHC
SIG-3523-26	Proliferating Cell Nuclear Antigen (PCNA) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	243	IHC

CatalogNo	Product Name	Size	EUR	Product Line
SIG-3610-1000	BRST-1 (CU-18) Monoclonal Antibody (SIGNET)	1.0 mL	688	IHC
SIG-3610-16	BRST-1 (CU-18) Monoclonal Antibody (SIGNET)	L1 6mL	445	IHC
SIG-3610-26	BRST-1 (CU-18) Monoclonal Antibody (SIGNET)	L2 6mL	416	IHC
SIG-3610-36	BRST-1 (CU-18) Monoclonal Antibody (SIGNET)	LV 6mL	458	IHC
SIG-3611-1000	GCDFP-15 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	688	IHC
SIG-3611-16	GCDFP-15 Monoclonal Antibody, Purified (SIGNET)	L1 6mL	445	IHC
SIG-3611-26	GCDFP-15 Monoclonal Antibody, Purified (SIGNET)	L2 6mL	416	IHC
SIG-3611-36	GCDFP-15 Monoclonal Antibody, Purified (SIGNET)	LV 6mL	458	IHC
SIG-3612-1000	BRST-3 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	688	IHC
SIG-3612-16	BRST-3 Monoclonal Antibody, Purified (SIGNET)	L1 6mL	445	IHC
SIG-3612-26	BRST-3 Monoclonal Antibody, Purified (SIGNET)	L2 6mL	416	IHC
SIG-3612-36	BRST-3 Monoclonal Antibody, Purified (SIGNET)	LV 6mL	458	IHC
SIG-3614-1000	CA 15-3 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	718	IHC
SIG-3615-26	CA 72.4 Monoclonal Antibody, Purified (SIGNET)	L2 6mL	342	IHC
SIG-3616-1000	CA 19-9 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	964	IHC
SIG-3616-16	CA 19-9 Monoclonal Antibody, Purified (SIGNET)	L1 6mL	458	IHC
SIG-3616-26	CA 19-9 Monoclonal Antibody, Purified (SIGNET)	L2 6mL	343	IHC
SIG-3617-1000	CA 125 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	646	IHC
SIG-3617-16	CA 125 Monoclonal Antibody, Purified (SIGNET)	L1 6mL	338	IHC
SIG-3617-26	CA 125 Monoclonal Antibody, Purified (SIGNET)	L2 6mL	301	IHC
SIG-3617-36	CA 125 Monoclonal Antibody, Purified (SIGNET)	LV 6mL	458	IHC
SIG-3619-1000	LK26 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	660	IHC
SIG-3620-1000	CD99 (Ewing's Sarcoma Marker) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	669	IHC
SIG-3620-16	CD99 (Ewing's Sarcoma Marker) Monoclonal Antibody, Purified (SIGNET)	L1 6mL	358	IHC
SIG-3620-26	CD99 (Ewing's Sarcoma Marker) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	315	IHC
SIG-3620-36	CD99 (Ewing's Sarcoma Marker) Monoclonal Antibody, Purified (SIGNET)	LV 6mL	458	IHC
SIG-3621-1000	MUC-1 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	688	IHC
SIG-3622-1000	Apo-D (GCDFP-24) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	688	IHC
SIG-3623-1000	CAK1 Monoclonal Antibody (SIGNET)	1.0 mL	811	IHC
SIG-3625-26	CD3 Polyclonal Antibody, Purified (SIGNET)	L2 6mL	214	IHC
SIG-3629-1000	CD8 Monoclonal Antibody (SIGNET)	1.0 mL	631	IHC
SIG-3629-26	CD8 Monoclonal Antibody (SIGNET)	L2 6mL	183	IHC
SIG-3630-1000	CD10/CALLA Monoclonal Antibody (SIGNET)	1.0 mL	791	IHC
SIG-3630-26	CD10/CALLA Monoclonal Antibody (SIGNET)	L2 6mL	271	IHC
SIG-3632-1000	CD31 (JC/70A) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	400	IHC
SIG-3632-26	CD31 (JC/70A) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	271	IHC
SIG-3633-1000	CD35 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	430	IHC
SIG-3633-26	CD35 Monoclonal Antibody, Purified (SIGNET)	L2 6mL	152	IHC
SIG-3634-1000	CD68 Monoclonal Antibody (SIGNET)	1.0 mL	372	IHC
SIG-3634-26	CD68 Monoclonal Antibody (SIGNET)	L2 6mL	185	IHC
SIG-3636-1000	CD23 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	372	IHC
SIG-3636-16	CD23 Monoclonal Antibody, Purified (SIGNET)	L1 6mL	185	IHC
SIG-3636-26	CD23 Monoclonal Antibody, Purified (SIGNET)	L2 6mL	128	IHC
SIG-3640-1000	CD43 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	395	IHC
SIG-3640-16	CD43 Monoclonal Antibody, Purified (SIGNET)	L1 6mL	229	IHC
SIG-3640-26	CD43 Monoclonal Antibody, Purified (SIGNET)	L2 6mL	185	IHC
SIG-3646-1000	CD45RB Monoclonal Antibody, Purified (SIGNET)	1.0 mL	343	IHC
SIG-3646-16	CD45RB Monoclonal Antibody, Purified (SIGNET)	L1 6mL	301	IHC
SIG-3646-26	CD45RB Monoclonal Antibody, Purified (SIGNET)	L2 6mL	271	IHC
SIG-3648-26	CD50 Monoclonal Antibody, Purified (SIGNET)	L2 6mL	315	IHC
SIG-3658-26	CD79 α Monoclonal Antibody (SIGNET)	L2 6mL	398	IHC
SIG-3660-1000	Estrogen Receptor (6F11) Monoclonal Antibody (SIGNET)	1.0 mL	1122	IHC
SIG-3665-1000	Estrogen Receptor (D75) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	548	IHC
SIG-3665-16	Estrogen Receptor (D75) Monoclonal Antibody, Purified (SIGNET)	L1 6mL	339	IHC
SIG-3665-26	Estrogen Receptor (D75) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	287	IHC
SIG-3670-1000	Estrogen Receptor (CO1531) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	548	IHC
SIG-3670-16	Estrogen Receptor (CO1531) Monoclonal Antibody, Purified (SIGNET)	L1 6mL	339	IHC
SIG-3670-26	Estrogen Receptor (CO1531) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	287	IHC
SIG-3701-1000	Cytokeratin 7 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	555	IHC
SIG-3701-26	Cytokeratin 7 Monoclonal Antibody, Purified (SIGNET)	L2 6mL	275	IHC
SIG-3710-26	Cytokeratin, LMW Monoclonal Antibody (SIGNET)	L2 6mL	217	IHC
SIG-3720-1000	Cytokeratin 20 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	1093	IHC
SIG-3720-26	Cytokeratin 20 Monoclonal Antibody, Purified (SIGNET)	L2 6mL	275	IHC
SIG-3725-26	Cytokeratin, HMW Monoclonal Antibody (SIGNET)	L2 6mL	217	IHC
SIG-3730-0100	D2-40 Lymphatic Endothelial Marker Monoclonal Antibody, Purified (SIGNET)	0.1 mL	407	IHC
SIG-3730-1000	D2-40 Lymphatic Endothelial Marker Monoclonal Antibody, Purified (SIGNET)	1.0 mL	714	IHC
SIG-3730-16	D2-40 Lymphatic Endothelial Marker Monoclonal Antibody, Purified (SIGNET)	L1 6mL	441	IHC
SIG-3730-26	D2-40 Lymphatic Endothelial Marker Monoclonal Antibody, Purified (SIGNET)	L2 6mL	379	IHC
SIG-3731-0100	HE4 Polyclonal Antibody (SIGNET)	0.1 mL	194	IHC
SIG-3731-1000	HE4 Polyclonal Antibody (SIGNET)	1.0 mL	683	IHC
SIG-3731-16	HE4 Polyclonal Antibody (SIGNET)	L1 6mL	413	IHC
SIG-3731-26	HE4 Polyclonal Antibody (SIGNET)	L2 6mL	348	IHC
SIG-3735-1000	HIP-1 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	741	IHC
SIG-3735-16	HIP-1 Monoclonal Antibody, Purified (SIGNET)	L1 6mL	458	IHC
SIG-3735-26	HIP-1 Monoclonal Antibody, Purified (SIGNET)	L2 6mL	395	IHC
SIG-3738-1000	Vascular Endothelial Growth Factor (VEGF) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	640	IHC
SIG-3738-16	Vascular Endothelial Growth Factor (VEGF) Monoclonal Antibody, Purified (SIGNET)	L1 6mL	486	IHC

CatalogNo	Product Name	Size	EUR	Product Line
SIG-3738-26	Vascular Endothelial Growth Factor (VEGF) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	456	IHC
SIG-3740-1000	Hepsin Polyclonal Antibody, Purified (SIGNET)	1.0 mL	741	IHC
SIG-3740-16	Hepsin Polyclonal Antibody, Purified (SIGNET)	L1 6mL	453	IHC
SIG-3740-26	Hepsin Polyclonal Antibody, Purified (SIGNET)	L2 6mL	380	IHC
SIG-3745-1000	Microphthalmia Transcription Factor (MITF) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	612	IHC
SIG-3745-16	Microphthalmia Transcription Factor (MITF) Monoclonal Antibody, Purified (SIGNET)	L1 6mL	379	IHC
SIG-3745-26	Microphthalmia Transcription Factor (MITF) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	325	IHC
SIG-3750-0100	CEACAM 6 Monoclonal Antibody (SIGNET)	0.1 mL	302	IHC
SIG-3750-0500	CEACAM 6 Monoclonal Antibody (SIGNET)	0.5 mL	1272	IHC
SIG-3751-0100	CEACAM 5 Monoclonal Antibody (SIGNET)	0.1 mL	302	IHC
SIG-3751-0500	CEACAM 5 Monoclonal Antibody (SIGNET)	0.5 mL	1272	IHC
SIG-3752-0100	CEACAM 5 Monoclonal Antibody (SIGNET)	0.1 mL	302	IHC
SIG-3752-0500	CEACAM 5 Monoclonal Antibody (SIGNET)	0.5 mL	1272	IHC
SIG-3753-0100	CEACAM 1,3,4,5,6 Monoclonal Antibody (SIGNET)	0.1 mL	302	IHC
SIG-3753-0500	CEACAM 1,3,4,5,6 Monoclonal Antibody (SIGNET)	0.5 mL	1263	IHC
SIG-3754-0100	CEACAM 5,6 Monoclonal Antibody (SIGNET)	0.1 mL	302	IHC
SIG-3754-0500	CEACAM 5,6 Monoclonal Antibody (SIGNET)	0.5 mL	1272	IHC
SIG-3755-0100	CEACAM 1,5,6,8 Monoclonal Antibody (SIGNET)	0.1 mL	302	IHC
SIG-3755-0500	CEACAM 1,5,6,8 Monoclonal Antibody (SIGNET)	0.5 mL	1272	IHC
SIG-3756-0100	CEACAM 1 Monoclonal Antibody (SIGNET)	0.1 mL	302	IHC
SIG-3756-0500	CEACAM 1 Monoclonal Antibody (SIGNET)	0.5 mL	1272	IHC
SIG-3760-1000	Bin-1 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	658	IHC
SIG-3760-0500	Bin-1 Monoclonal Antibody, Purified (SIGNET)	0.5 mL	374	IHC
SIG-3770-1000	Endoglin (CD105) Clone SN6h Monoclonal Antibody, Purified (Signet)	1.0 mL	600	IHC
SIG-3770-26	Endoglin (CD105) Clone SN6h Monoclonal Antibody, Purified (Signet)	L2 6mL	392	IHC
SIG-38001-kit	OPD Diluent and Tablets	kit	93	Dedham Neuro
SIG-38010-1000	URO-1 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	448	IHC
SIG-3801-26	Androgen Receptor Monoclonal Antibody, Purified (SIGNET)	L2 6mL	275	IHC
SIG-38020-1000	URO-2 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	461	IHC
SIG-38040-1000	URO-4 (DPP-IV, CD26, ABP-26) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	461	IHC
SIG-38050-1000	URO-5 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	461	IHC
SIG-38070-1000	URO-7 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	461	IHC
SIG-38080-1000	URO-8 Monoclonal Antibody (SIGNET)	1.0 mL	461	IHC
SIG-38090-1000	URO-9 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	461	IHC
SIG-38110-1000	MEL-1 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	439	IHC
SIG-38130-1000	MEL-3 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	439	IHC
SIG-38140-1000	MEL-4 Monoclonal Antibody (SIGNET)	1.0 mL	439	IHC
SIG-38150-1000	MEL-5 Monoclonal Antibody (SIGNET)	1.0 mL	439	IHC
SIG-38150-26	MEL-5 Monoclonal Antibody (SIGNET)	L2 6mL	275	IHC
SIG-38150-36	MEL-5 Monoclonal Antibody (SIGNET)	LV 6mL	398	IHC
SIG-38160-1000	MART-1 (M2-7C10) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	654	IHC
SIG-38160-26	MART-1 (M2-7C10) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	482	IHC
SIG-38160-36	MART-1 (M2-7C10) Monoclonal Antibody, Purified (SIGNET)	LV 6mL	482	IHC
SIG-38165-1000	MART-1 (M2-9E3) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	654	IHC
SIG-3820-26	TdT Polyclonal Antibody, Purified (SIGNET)	L2 6mL	349	IHC
SIG-3830-26	Thyroid Transcription Factor (TTF-1) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	349	IHC
SIG-3840-1000	Tubulin (TU-20) Monoclonal Antibody (SIGNET)	1.0 mL	749	IHC
SIG-3840-26	Tubulin (TU-20) Monoclonal Antibody (SIGNET)	L2 6mL	459	IHC
SIG-3860-1000	Epidermal Growth Factor Receptor (111.6) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	820	IHC
SIG-3860-26	Epidermal Growth Factor Receptor (111.6) Monoclonal Antibody, Purified (SIGNET)	L2 6mL	459	IHC
SIG-38710-1000	C219 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	724	IHC
SIG-38720-1000	C494 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	838	IHC
SIG-38730-1000	JSB1 Monoclonal Antibody (SIGNET)	1.0 mL	586	IHC
SIG-38740-1000	4E3 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	838	IHC
SIG-38750-1000	Lung Related Protein Monoclonal Antibody (SIGNET)	1.0 mL	691	IHC
SIG-38760-1000	Multidrug Resistance Associated Protein (MRPr1) Monoclonal Antibody (SIGNET)	1.0 mL	848	IHC
SIG-38765-1000	Multidrug Resistance Protein (QCRL1) Monoclonal Antibody (SIGNET)	1.0 mL	838	IHC
SIG-38767-1000	Multidrug Resistance Protein (QCRL3) Monoclonal Antibody (SIGNET)	1.0 mL	838	IHC
SIG-38770-1000	Multidrug Resistance Protein (MRPm6) Monoclonal Antibody (SIGNET)	1.0 mL	765	IHC
SIG-38775-1000	Multidrug Resistance Protein 1 (MRP1) Monoclonal Antibody (SIGNET)	1.0 mL	765	IHC
SIG-38780-1000	cMOAT/MRP2 (M2I-4) Monoclonal Antibody (SIGNET)	1.0 mL	765	IHC
SIG-38785-1000	cMOAT/MRP2 (M2III-6) Monoclonal Antibody (SIGNET)	1.0 mL	765	IHC
SIG-38790-1000	Lung Resistance Related Protein (LRP) Monoclonal Antibody (SIGNET)	1.0 mL	765	IHC
SIG-38795-1000	Multidrug Resistance Protein 3 (M3II-9) Monoclonal Antibody (SIGNET)	1.0 mL	765	IHC
SIG-38800-1000	Multidrug Resistance Protein 3 (M3II-21) Monoclonal Antibody (SIGNET)	1.0 mL	765	IHC
SIG-38805-1000	cMOAT/MRP2 (M2II-12) Monoclonal Antibody (SIGNET)	1.0 mL	765	IHC
SIG-38810-1000	Multidrug Resistance Protein 5 (M5I-1) Monoclonal Antibody (SIGNET)	1.0 mL	765	IHC
SIG-38815-1000	Multidrug Resistance Protein 5 (M5II-54) Monoclonal Antibody (SIGNET)	1.0 mL	765	IHC
SIG-38820-1000	Breast Cancer Resistance Protein (BXP-34) Monoclonal Antibody (SIGNET)	1.0 mL	765	IHC
SIG-38822-1000	Breast Cancer Resistance Protein (BXP-21) Monoclonal Antibody (SIGNET)	1.0 mL	765	IHC
SIG-38825-1000	Multidrug Resistance Protein 6 (M6II-7) Monoclonal Antibody (SIGNET)	1.0 mL	765	IHC
SIG-38827-1000	Multidrug Resistance Protein 6 (M6II-21) Monoclonal Antibody (SIGNET)	1.0 mL	765	IHC
SIG-38830-1000	Multidrug Resistance Protein 6 (M6II-31) Monoclonal Antibody (SIGNET)	1.0 mL	765	IHC
SIG-38910-pair	Apolipoprotein-D (ApoD) Matched Pair Monoclonal Antibodies (SIGNET)	pair	2760	Dedham Neuro
SIG-38930-050	Human Apolipoprotein D Protein (SIGNET)	50 uL	737	Dedham Neuro
SIG-38940-kit	Beta-Amyloid 1-40 ELISA Kit (SIGNET)	kit	692	Dedham Neuro
SIG-38942-kit	Beta-Amyloid 1-42 ELISA Kit (SIGNET)	kit	692	Dedham Neuro

CatalogNo	Product Name	Size	EUR	Product Line
SIG-38950-kit	BetaMark™ Beta-Amyloid x-40 ELISA Kit	kit	764	Dedham Neuro
SIG-38952-kit	BetaMark™ Beta-Amyloid x-42 ELISA Kit	kit	764	Dedham Neuro
SIG-38954-kit	Colorimetric BetaMark Beta-Amyloid x-40 ELISA Kit	kit	732	Dedham Neuro
SIG-38956-kit	Colorimetric BetaMark Beta-Amyloid x-42 ELISA Kit	kit	732	Dedham Neuro
SIG-38960-kit	sAPP Beta ELISA kit	kit	795	Added April 10th
SIG-3903-1000	Cyclin E Monoclonal Antibody, Purified (SIGNET)	1.0 mL	612	IHC
SIG-3903-16	Cyclin E Monoclonal Antibody, Purified (SIGNET)	L1 6mL	379	IHC
SIG-3903-26	Cyclin E Monoclonal Antibody, Purified (SIGNET)	L2 6mL	325	IHC
SIG-3905-1000	Cyclin D1 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	640	IHC
SIG-3905-16	Cyclin D1 Monoclonal Antibody, Purified (SIGNET)	L1 6mL	395	IHC
SIG-3905-26	Cyclin D1 Monoclonal Antibody, Purified (SIGNET)	L2 6mL	333	IHC
SIG-3906-1000	Cyclin D3 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	589	IHC
SIG-39061-050	NMDA Receptor NR2A Polyclonal Antibody (SIGNET)	50 ug	479	Dedham Neuro
SIG-3906-16	Cyclin D3 Monoclonal Antibody, Purified (SIGNET)	L1 6mL	365	IHC
SIG-3906-26	Cyclin D3 Monoclonal Antibody, Purified (SIGNET)	L2 6mL	313	IHC
SIG-39063-050	NMDA Receptor NR2B Polyclonal Antibody (SIGNET)	50 ug	479	Dedham Neuro
SIG-39070-1000	MLK-2 Polyclonal Antibody (SIGNET)	1.0 mL	408	Dedham Neuro
SIG-39080-050	Apo-J/Clusterin Polyclonal Antibody (SIGNET)	50 ug	581	Dedham Neuro
SIG-39085-050	Apo-J Protein, Purified	50 uL	720	Dedham Neuro
SIG-39110-1000	β-Amyloid (1-8) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	1828	Dedham Neuro
SIG-39110-0200	β-Amyloid (1-8) Monoclonal Antibody, Purified (SIGNET)	0.2 mL	581	Dedham Neuro
SIG-39110-0500	β-Amyloid (1-8) Monoclonal Antibody, Purified (SIGNET)	0.5 mL	1115	Dedham Neuro
SIG-3911-1000	L1 (CD171) Monoclonal Antibody (SIGNET)	1.0 mL	683	IHC
SIG-3911-16	L1 (CD171) Monoclonal Antibody (SIGNET)	L1 6mL	413	IHC
SIG-3911-26	L1 (CD171) Monoclonal Antibody (SIGNET)	L2 6mL	348	IHC
SIG-39120-1000	β-Amyloid (1-8) Monoclonal Antibody (SIGNET)	1.0 mL	1828	Dedham Neuro
SIG-39120-0200	β-Amyloid (1-8) Monoclonal Antibody (SIGNET)	0.2 mL	581	Dedham Neuro
SIG-39120-0500	β-Amyloid (1-8) Monoclonal Antibody (SIGNET)	0.5 mL	1115	Dedham Neuro
SIG-3912-1000	Phosphatase and Tensin Homolog (PTEN) Monoclonal Antibody (SIGNET)	1.0 mL	683	IHC
SIG-3912-16	Phosphatase and Tensin Homolog (PTEN) Monoclonal Antibody (SIGNET)	L1 6mL	413	IHC
SIG-3912-26	Phosphatase and Tensin Homolog (PTEN) Monoclonal Antibody (SIGNET)	L2 6mL	348	IHC
SIG-39130-050	β-Amyloid 40 Polyclonal Antibody, Purified (SIGNET)	50 ug	741	Dedham Neuro
SIG-39131-1000	β-Amyloid 40 Polyclonal Antibody, Purified IgG Fraction (SIGNET)	1.0 mL	1828	Dedham Neuro
SIG-39131-0200	β-Amyloid 40 Polyclonal Antibody, Purified IgG Fraction (SIGNET)	0.2 mL	581	Dedham Neuro
SIG-39131-0500	β-Amyloid 40 Polyclonal Antibody, Purified IgG Fraction (SIGNET)	0.5 mL	1115	Dedham Neuro
SIG-39132-1000	β-Amyloid 40 Polyclonal Antibody, Biotin Labeled IgG Fraction (SIGNET)	1.0 mL	1384	Dedham Neuro
SIG-39132-0200	β-Amyloid 40 Polyclonal Antibody, Biotin Labeled IgG Fraction (SIGNET)	0.2 mL	714	Dedham Neuro
SIG-39132-0500	β-Amyloid 40 Polyclonal Antibody, Biotin Labeled IgG Fraction (SIGNET)	0.5 mL	790	Dedham Neuro
SIG-39133-050	β-Amyloid 40 Polyclonal Antibody, Biotin Labeled (SIGNET)	50 ug	790	Dedham Neuro
SIG-39134-1000	β-Amyloid 42 Polyclonal Antibody, Purified IgG Fraction (SIGNET)	1.0 mL	1828	Dedham Neuro
SIG-39134-0200	β-Amyloid 42 Polyclonal Antibody, Purified IgG Fraction (SIGNET)	0.2 mL	581	Dedham Neuro
SIG-39134-0500	β-Amyloid 42 Polyclonal Antibody, Purified IgG Fraction (SIGNET)	0.5 mL	1115	Dedham Neuro
SIG-39135-1000	β-Amyloid 42 Polyclonal Antibody, Biotin Labeled IgG Fraction (SIGNET)	1.0 mL	2427	Dedham Neuro
SIG-39135-0200	β-Amyloid 42 Polyclonal Antibody, Biotin Labeled IgG Fraction (SIGNET)	0.2 mL	714	Dedham Neuro
SIG-39135-0500	β-Amyloid 42 Polyclonal Antibody, Biotin Labeled IgG Fraction (SIGNET)	0.5 mL	1384	Dedham Neuro
SIG-39136-050	β-Amyloid 42 Polyclonal Antibody, Purified (SIGNET)	50 ug	741	Dedham Neuro
SIG-39137-050	β-Amyloid 42 Polyclonal Antibody, Biotin Labeled (SIGNET)	50 ug	790	Dedham Neuro
SIG-39138-050	sAPPβ Polyclonal Antibody, Purified	0.05 mg	714	Dedham Neuro
SIG-39139-050	sAPPα Polyclonal Antibody, Purified	0.05 mg	714	Dedham Neuro
SIG-39140-1000	β-Amyloid 1-40 (11A50-B10) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	1828	Dedham Neuro
SIG-39140-0200	β-Amyloid 1-40 (11A50-B10) Monoclonal Antibody, Purified (SIGNET)	0.2 mL	581	Dedham Neuro
SIG-39140-0500	β-Amyloid 1-40 (11A50-B10) Monoclonal Antibody, Purified (SIGNET)	0.5 mL	1115	Dedham Neuro
SIG-39141-1000	β-Amyloid 1-40 (11A50-B10) Monoclonal Antibody, Biotin Labeled (SIGNET)	1.0 mL	2209	Dedham Neuro
SIG-39141-0200	β-Amyloid 1-40 (11A50-B10) Monoclonal Antibody, Biotin Labeled (SIGNET)	0.2 mL	651	Dedham Neuro
SIG-39141-0500	β-Amyloid 1-40 (11A50-B10) Monoclonal Antibody, Biotin Labeled (SIGNET)	0.5 mL	1261	Dedham Neuro
SIG-39142-1000	β-Amyloid 1-42 (12F4) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	1828	Dedham Neuro
SIG-39142-0200	β-Amyloid 1-42 (12F4) Monoclonal Antibody, Purified (SIGNET)	0.2 mL	581	Dedham Neuro
SIG-39142-0500	β-Amyloid 1-42 (12F4) Monoclonal Antibody, Purified (SIGNET)	0.5 mL	1113	Dedham Neuro
SIG-39143-1000	β-Amyloid 1-43 (9C4) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	1828	Dedham Neuro
SIG-39143-0200	β-Amyloid 1-43 (9C4) Monoclonal Antibody, Purified (SIGNET)	0.2 mL	581	Dedham Neuro
SIG-39143-0500	β-Amyloid 1-43 (9C4) Monoclonal Antibody, Purified (SIGNET)	0.5 mL	1115	Dedham Neuro
SIG-39144-1000	β-Amyloid 1-42 (12F4) Monoclonal Antibody, Biotin Labeled (SIGNET)	1.0 mL	2209	Dedham Neuro
SIG-39144-0200	β-Amyloid 1-42 (12F4) Monoclonal Antibody, Biotin Labeled (SIGNET)	0.2 mL	651	Dedham Neuro
SIG-39144-0500	β-Amyloid 1-42 (12F4) Monoclonal Antibody, Biotin Labeled (SIGNET)	0.5 mL	1261	Dedham Neuro
SIG-39145-1000	β-Amyloid 1-43 (9C4) Monoclonal Antibody, Biotin Labeled (SIGNET)	1.0 mL	2209	Dedham Neuro
SIG-39145-0200	β-Amyloid 1-43 (9C4) Monoclonal Antibody, Biotin Labeled (SIGNET)	0.2 mL	651	Dedham Neuro
SIG-39145-0500	β-Amyloid 1-43 (9C4) Monoclonal Antibody, Biotin Labeled (SIGNET)	0.5 mL	1261	Dedham Neuro
SIG-39146-0200	Beta Amyloid 1-40 (11A50-B10) Monoclonal Antibody, HRP Labeled (SIGNET)	0.2 mL	692	Dedham Neuro
SIG-39147-0200	β-Amyloid 1-42 (12F4) Monoclonal Antibody, HRP Labeled (SIGNET)	0.2 mL	692	Dedham Neuro
SIG-39148-050	AICD (APP Intracellular Domain) Polyclonal Antibody	50 ug	581	Dedham Neuro
SIG-39150-050	Amyloid Precursor Protein (APP) C-Terminal Fragment Polyclonal Antibody, Purified (SIGNET)	50 ug	581	Dedham Neuro
SIG-3915-1000	ICBP90 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	603	IHC
SIG-39151-1000	β-Amyloid, Rodent Specific Polyclonal Antibody, IgG Fraction (SIGNET)	1.0 mL	2427	Dedham Neuro
SIG-39151-0200	β-Amyloid, Rodent Specific Polyclonal Antibody, IgG Fraction (SIGNET)	0.2 mL	714	Dedham Neuro
SIG-39151-0500	β-Amyloid, Rodent Specific Polyclonal Antibody, IgG Fraction (SIGNET)	0.5 mL	1384	Dedham Neuro
SIG-39153-050	β-Amyloid, Rodent Specific Polyclonal Antibody, Purified (SIGNET)	50 ug	714	Dedham Neuro
SIG-39154-050	β-Amyloid, Rodent Specific Polyclonal Antibody, Biotin Labeled (SIGNET)	50 ug	754	Dedham Neuro

CatalogNo	Product Name	Size	EUR	Product Line
SIG-3915-0500	ICBP90 Monoclonal Antibody, Purified (SIGNET)	0.5 mL	374	IHC
SIG-39160-050	Beta Amyloid 38 Clone BA1-13 Monoclonal Antibody, Purified	50ug	480	Dedham Neuro
SIG-39164-050	Beta Amyloid 40 Rabbit Monoclonal, Purified	50ug	480	Added April 10th
SIG-39168-050	Beta Amyloid 42 Clone BA3-9 Monoclonal Antibody, Purified	50ug	480	Dedham Neuro
SIG-39172-0100	Notch Polyclonal Antibody	100ug	559	Dedham Neuro
SIG-39180-1000	Amyloid Precursor Protein (APP) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	1825	Dedham Neuro
SIG-39180-0200	Amyloid Precursor Protein (APP) Monoclonal Antibody, Purified (SIGNET)	0.2 mL	587	Dedham Neuro
SIG-39180-0500	Amyloid Precursor Protein (APP) Monoclonal Antibody, Purified (SIGNET)	0.5 mL	1163	Dedham Neuro
SIG-39181-0200	Amyloid Precursor Protein (APP) Monoclonal Antibody, Biotinylated	0.2 mL	687	Dedham Neuro
SIG-39184-1000	Amyloid Precursor protein (APP) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	1825	Dedham Neuro
SIG-39184-0200	Amyloid Precursor protein (APP) Monoclonal Antibody, Purified (SIGNET)	0.2 mL	587	Dedham Neuro
SIG-39184-0500	Amyloid Precursor protein (APP) Monoclonal Antibody, Purified (SIGNET)	0.5 mL	1163	Dedham Neuro
SIG-39188-1000	Amyloid Precursor Protein (APP) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	1825	Dedham Neuro
SIG-39188-0200	Amyloid Precursor Protein (APP) Monoclonal Antibody, Purified (SIGNET)	0.2 mL	587	Dedham Neuro
SIG-39188-0500	Amyloid Precursor Protein (APP) Monoclonal Antibody, Purified (SIGNET)	0.5 mL	1163	Dedham Neuro
SIG-39190-1000	Presenilin 1 (NT) Monoclonal Antibody, Purified	1.0mL	1757	Dedham Neuro
SIG-39190-0200	Presenilin 1 (NT) Monoclonal Antibody, Purified	0.2mL	554	Dedham Neuro
SIG-39190-0500	Presenilin 1 (NT) Monoclonal Antibody, Purified	0.5mL	1120	Dedham Neuro
SIG-39194-1000	Presenilin 1 (NT1) Monoclonal Antibody, Purified	1.0 mL	1757	Dedham Neuro
SIG-39194-0200	Presenilin 1 (NT1) Monoclonal Antibody, Purified	0.2 mL	554	Dedham Neuro
SIG-39194-0500	Presenilin 1 (NT1) Monoclonal Antibody, Purified	0.5 mL	1120	Dedham Neuro
SIG-39198-1000	Presenilin 1 LOOP Monoclonal Antibody, Purified	1.0 mL	1757	Dedham Neuro
SIG-39198-0200	Presenilin 1 LOOP Monoclonal Antibody, Purified	0.2 mL	554	Dedham Neuro
SIG-39198-0500	Presenilin 1 LOOP Monoclonal Antibody, Purified	0.5 mL	1120	Dedham Neuro
SIG-39200-1000	β-Amyloid, 17-24 (4G8) Monoclonal Antibody (SIGNET) <I>Exclusively Provided b	1.0 mL	3800	Dedham Neuro
SIG-39200-0200	β-Amyloid, 17-24 (4G8) Monoclonal Antibody (SIGNET) <I>Exclusively Provided b	0.2 mL	1134	Dedham Neuro
SIG-39200-0500	β-Amyloid, 17-24 (4G8) Monoclonal Antibody (SIGNET) <I>Exclusively Provided b	0.5 mL	2278	Dedham Neuro
SIG-39220-1000	β-Amyloid, 17-24 (4G8) Monoclonal Antibody, Purified (SIGNET) <I>Exclusively P	1.0 mL	1897	Dedham Neuro
SIG-39220-0200	β-Amyloid, 17-24 (4G8) Monoclonal Antibody, Purified (SIGNET) <I>Exclusively P	0.2 mL	609	Dedham Neuro
SIG-39220-0500	β-Amyloid, 17-24 (4G8) Monoclonal Antibody, Purified (SIGNET) <I>Exclusively P	0.5 mL	1163	Dedham Neuro
SIG-39240-1000	β-Amyloid, 17-24 (4G8) Monoclonal Antibody, Biotin Labeled (SIGNET) <I>Exclus	1.0 mL	2427	Dedham Neuro
SIG-39240-0200	β-Amyloid, 17-24 (4G8) Monoclonal Antibody, Biotin Labeled (SIGNET) <I>Exclus	0.2 mL	714	Dedham Neuro
SIG-39240-0500	β-Amyloid, 17-24 (4G8) Monoclonal Antibody, Biotin Labeled (SIGNET) <I>Exclus	0.5 mL	1384	Dedham Neuro
SIG-39245-0200	β-Amyloid, 17-24 (4G8) Monoclonal Antibody, HRP Labeled (SIGNET) <I>Exclusi	0.2 mL	719	Dedham Neuro
SIG-39300-1000	β-Amyloid, 1-16 (6E10) Monoclonal Antibody (SIGNET) <I>Exclusively Provided b	1.0 mL	3800	Dedham Neuro
SIG-39300-0200	β-Amyloid, 1-16 (6E10) Monoclonal Antibody (SIGNET) <I>Exclusively Provided b	0.2 mL	1134	Dedham Neuro
SIG-39300-0500	β-Amyloid, 1-16 (6E10) Monoclonal Antibody (SIGNET) <I>Exclusively Provided b	0.5 mL	2278	Dedham Neuro
SIG-39320-1000	β-Amyloid, 1-16 (6E10) Monoclonal Antibody, Purified (SIGNET) <I>Exclusively P	1.0 mL	1897	Dedham Neuro
SIG-39320-0200	β-Amyloid, 1-16 (6E10) Monoclonal Antibody, Purified (SIGNET) <I>Exclusively P	0.2 mL	609	Dedham Neuro
SIG-39320-0500	β-Amyloid, 1-16 (6E10) Monoclonal Antibody, Purified (SIGNET) <I>Exclusively P	0.5 mL	1163	Dedham Neuro
SIG-39321-1000	Beta Amyloid, 1-16 (6E10) Monoclonal Antibody Low Endotoxin, Purified	1.0 mL	2271	
SIG-39322-050	β-Amyloid, 1-16 Polyclonal Antibody, Purified	50 ug	559	Dedham Neuro
SIG-39330-0200	β-Amyloid, 1-16 (6E10) Monoclonal Antibody, F(Ab)2 Fragments (SIGNET) <I>Ex	0.2 mL	719	Dedham Neuro
SIG-39340-1000	β-Amyloid, 1-16 (6E10) Monoclonal Antibody, Biotin Labeled (SIGNET) <I>Exclusi	1.0 mL	2427	Dedham Neuro
SIG-39340-0200	β-Amyloid, 1-16 (6E10) Monoclonal Antibody, Biotin Labeled (SIGNET) <I>Exclusi	0.2 mL	714	Dedham Neuro
SIG-39340-0500	β-Amyloid, 1-16 (6E10) Monoclonal Antibody, Biotin Labeled (SIGNET) <I>Exclusi	0.5 mL	1384	Dedham Neuro
SIG-39345-0200	β-Amyloid, 1-16 (6E10) Monoclonal Antibody, HRP Labeled (SIGNET) <I>Exclusi	0.2 mL	719	Dedham Neuro
SIG-39347-0200	β-Amyloid, 1-16 (6E10) Monoclonal Antibody, Alexa Fluor® 488 Labeled (SIG	0.2 mL	740	Dedham Neuro
SIG-39350-1000	β-Amyloid (20.1) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	1825	Dedham Neuro
SIG-39350-0200	β-Amyloid (20.1) Monoclonal Antibody, Purified (SIGNET)	0.2 mL	587	Dedham Neuro
SIG-39360-1000	Presenilin 1 Polyclonal Antibody (SIGNET)	1.0 mL	2007	Dedham Neuro
SIG-39360-0200	Presenilin 1 Polyclonal Antibody (SIGNET)	0.2 mL	636	Dedham Neuro
SIG-39360-0500	Presenilin 1 Polyclonal Antibody (SIGNET)	0.5 mL	1225	Dedham Neuro
SIG-39370-0200	Nicastrin Polyclonal Antibody, Purified (SIGNET)	0.2 mL	636	Dedham Neuro
SIG-39400-1000	Ubiquitin (5-25) Monoclonal Antibody (SIGNET)	1.0 mL	4027	Dedham Neuro
SIG-39400-0200	Ubiquitin (5-25) Monoclonal Antibody (SIGNET)	0.2 mL	1225	Dedham Neuro
SIG-39400-0500	Ubiquitin (5-25) Monoclonal Antibody (SIGNET)	0.5 mL	2427	Dedham Neuro
SIG-39410-1000	Tau (C3) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	1757	Dedham Neuro
SIG-39410-0200	Tau (C3) Monoclonal Antibody, Purified (SIGNET)	0.2 mL	454	Dedham Neuro
SIG-39410-0500	Tau (C3) Monoclonal Antibody, Purified (SIGNET)	0.5 mL	947	Dedham Neuro
SIG-39413-1000	Tau (Tau 5) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	1757	Dedham Neuro
SIG-39413-0200	Tau (Tau 5) Monoclonal Antibody, Purified (SIGNET)	0.2 mL	454	Dedham Neuro
SIG-39413-0500	Tau (Tau 5) Monoclonal Antibody, Purified (SIGNET)	0.5 mL	947	Dedham Neuro
SIG-39416-1000	Tau (Tau 12) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	1757	Dedham Neuro
SIG-39416-0200	Tau (Tau 12) Monoclonal Antibody, Purified (SIGNET)	0.2 mL	454	Dedham Neuro
SIG-39416-0500	Tau (Tau 12) Monoclonal Antibody, Purified (SIGNET)	0.5 mL	947	Dedham Neuro
SIG-39423-1000	Tau (Tau46) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	1559	Dedham Neuro
SIG-39423-0200	Tau (Tau46) Monoclonal Antibody, Purified (SIGNET)	0.2 mL	421	Dedham Neuro
SIG-39423-0500	Tau (Tau46) Monoclonal Antibody, Purified (SIGNET)	0.5 mL	893	Dedham Neuro
SIG-39426-1000	Tau (Tau 2) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	1757	Dedham Neuro
SIG-39426-0200	Tau (Tau 2) Monoclonal Antibody, Purified (SIGNET)	0.2 mL	454	Dedham Neuro
SIG-39426-0500	Tau (Tau 2) Monoclonal Antibody, Purified (SIGNET)	0.5 mL	947	Dedham Neuro
SIG-39430-1000	Paired Helical Filaments (PHF 6) Monoclonal Antibody, Purified	1.0 mL	1559	Dedham Neuro
SIG-39430-0200	Paired Helical Filaments (PHF 6) Monoclonal Antibody, Purified	0.2 mL	421	Dedham Neuro
SIG-39430-0500	Paired Helical Filaments (PHF 6) Monoclonal Antibody, Purified	0.5 mL	893	Dedham Neuro
SIG-39433-1000	Paired Helical Filaments (PHF 13) Monoclonal Antibody, Purified	1.0 mL	1620	Dedham Neuro
SIG-39433-0200	Paired Helical Filaments (PHF 13) Monoclonal Antibody, Purified	0.2 mL	436	Dedham Neuro

CatalogNo	Product Name	Size	EUR	Product Line
SIG-39433-0500	Paired Helical Filaments (PHF 13) Monoclonal Antibody, Purified	0.5 mL	928	Dedham Neuro
SIG-39436-1000	Nitrated Tau (Tau-nY18) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	1757	Dedham Neuro
SIG-39436-0200	Nitrated Tau (Tau-nY18) Monoclonal Antibody, Purified (SIGNET)	0.2 mL	454	Dedham Neuro
SIG-39436-0500	Nitrated Tau (Tau-nY18) Monoclonal Antibody, Purified (SIGNET)	0.5 mL	947	Dedham Neuro
SIG-39439-1000	Nitrated Tau (Tau-nY29) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	1757	Dedham Neuro
SIG-39439-0200	Nitrated Tau (Tau-nY29) Monoclonal Antibody, Purified (SIGNET)	0.2 mL	454	Dedham Neuro
SIG-39439-0500	Nitrated Tau (Tau-nY29) Monoclonal Antibody, Purified (SIGNET)	0.5 mL	947	Dedham Neuro
SIG-39500-1000	Ubiquitin (3-39) Monoclonal Antibody (SIGNET)	1.0 mL	4027	Dedham Neuro
SIG-39500-0200	Ubiquitin (3-39) Monoclonal Antibody (SIGNET)	0.2 mL	1225	Dedham Neuro
SIG-39500-0500	Ubiquitin (3-39) Monoclonal Antibody (SIGNET)	0.5 mL	2427	Dedham Neuro
SIG-39523-1000	Parkin (Prk 109) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	1620	Dedham Neuro
SIG-39523-0200	Parkin (Prk 109) Monoclonal Antibody, Purified (SIGNET)	0.2 mL	436	Dedham Neuro
SIG-39523-0500	Parkin (Prk 109) Monoclonal Antibody, Purified (SIGNET)	0.5 mL	928	Dedham Neuro
SIG-39530-1000	Parkin (Prk 8) Monoclonal Antibody, Purified	1.0 mL	1620	Dedham Neuro
SIG-39530-0200	Parkin (Prk 8) Monoclonal Antibody, Purified	0.2 mL	436	Dedham Neuro
SIG-39530-0500	Parkin (Prk 8) Monoclonal Antibody, Purified	0.5 mL	928	Dedham Neuro
SIG-39600-1000	Prion (3F4) Monoclonal Antibody (SIGNET) <I>Exclusively Provided by Covance</I>	1.0 mL	3800	Dedham Neuro
SIG-39600-0200	Prion (3F4) Monoclonal Antibody (SIGNET) <I>Exclusively Provided by Covance</I>	0.2 mL	1134	Dedham Neuro
SIG-39600-0500	Prion (3F4) Monoclonal Antibody (SIGNET) <I>Exclusively Provided by Covance</I>	0.5 mL	2278	Dedham Neuro
SIG-39620-1000	Prion (3F4) Monoclonal Antibody, Purified (SIGNET) <I>Exclusively Provided by Covar	1.0 mg	1897	Dedham Neuro
SIG-39620-0200	Prion (3F4) Monoclonal Antibody, Purified (SIGNET) <I>Exclusively Provided by Covar	0.2 mg	609	Dedham Neuro
SIG-39620-0500	Prion (3F4) Monoclonal Antibody, Purified (SIGNET) <I>Exclusively Provided by Covar	0.5 mg	1163	Dedham Neuro
SIG-39640-1000	Prion (3F4) Monoclonal Antibody, Biotin Labeled (SIGNET) <I>Exclusively Provided by	1.0 mg	2427	Dedham Neuro
SIG-39640-0200	Prion (3F4) Monoclonal Antibody, Biotin Labeled (SIGNET) <I>Exclusively Provided by	0.2 mg	714	Dedham Neuro
SIG-39640-0500	Prion (3F4) Monoclonal Antibody, Biotin Labeled (SIGNET) <I>Exclusively Provided by	0.5 mg	1384	Dedham Neuro
SIG-39700-1000	Alpha-I Spectrin Monoclonal Antibody (SIGNET)	1.0 mL	2007	Dedham Neuro
SIG-39700-0200	Alpha-I Spectrin Monoclonal Antibody (SIGNET)	0.2 mL	636	Dedham Neuro
SIG-39700-0500	Alpha-I Spectrin Monoclonal Antibody (SIGNET)	0.5 mL	1225	Dedham Neuro
SIG-39702-1000	Alpha-II Spectrin Monoclonal Antibody (SIGNET)	1.0 mL	2007	Dedham Neuro
SIG-39702-0200	Alpha-II Spectrin Monoclonal Antibody (SIGNET)	0.2 mL	636	Dedham Neuro
SIG-39702-0500	Alpha-II Spectrin Monoclonal Antibody (SIGNET)	0.5 mL	1225	Dedham Neuro
SIG-39710-1000	Spectrin Binding Protein Monoclonal Antibody (SIGNET)	1.0 mL	2007	Dedham Neuro
SIG-39710-0200	Spectrin Binding Protein Monoclonal Antibody (SIGNET)	0.2 mL	636	Dedham Neuro
SIG-39710-0500	Spectrin Binding Protein Monoclonal Antibody (SIGNET)	0.5 mL	1225	Dedham Neuro
SIG-39720-1000	α-Synuclein (4D6) Monoclonal Antibody (SIGNET)	1.0 mL	1828	Dedham Neuro
SIG-39720-0200	α-Synuclein (4D6) Monoclonal Antibody (SIGNET)	0.2 mL	581	Dedham Neuro
SIG-39720-0500	α-Synuclein (4D6) Monoclonal Antibody (SIGNET)	0.5 mL	1115	Dedham Neuro
SIG-39725-1000	α-Synuclein (LB509) Monoclonal Antibody (SIGNET)	1.0 mL	1828	Dedham Neuro
SIG-39725-0200	α-Synuclein (LB509) Monoclonal Antibody (SIGNET)	0.2 mL	581	Dedham Neuro
SIG-39725-0500	α-Synuclein (LB509) Monoclonal Antibody (SIGNET)	0.5 mL	1115	Dedham Neuro
SIG-39726-1000	α-Synuclein (LB509) Monoclonal Antibody	1.0mL	2098	Dedham Neuro
SIG-39726-0200	α-Synuclein (LB509) Monoclonal Antibody	0.2mL	664	Dedham Neuro
SIG-39726-0500	α-Synuclein (LB509) Monoclonal Antibody	0.5mL	1279	Dedham Neuro
SIG-39730-1000	α-Synuclein (4B12) Monoclonal Antibody (SIGNET)	1.0 mL	1828	Dedham Neuro
SIG-39730-0200	α-Synuclein (4B12) Monoclonal Antibody (SIGNET)	0.2 mL	581	Dedham Neuro
SIG-39730-0500	α-Synuclein (4B12) Monoclonal Antibody (SIGNET)	0.5 mL	1115	Dedham Neuro
SIG-39740-1000	Apolipoprotein-E (D6E10) Monoclonal Antibody (SIGNET)	1.0 mL	1862	Dedham Neuro
SIG-39740-0200	Apolipoprotein-E (D6E10) Monoclonal Antibody (SIGNET)	0.2 mL	617	Dedham Neuro
SIG-39740-0500	Apolipoprotein-E (D6E10) Monoclonal Antibody (SIGNET)	0.5 mL	1134	Dedham Neuro
SIG-39750-1000	Apolipoprotein-E (E6D7) Monoclonal Antibody (SIGNET)	1.0 mL	1862	Dedham Neuro
SIG-39750-0200	Apolipoprotein-E (E6D7) Monoclonal Antibody (SIGNET)	0.2 mL	617	Dedham Neuro
SIG-39750-0500	Apolipoprotein-E (E6D7) Monoclonal Antibody (SIGNET)	0.5 mL	1134	Dedham Neuro
SIG-39760-1000	ERAB (10D12) Monoclonal Antibody (SIGNET)	1.0 mL	2007	Dedham Neuro
SIG-39760-0200	ERAB (10D12) Monoclonal Antibody (SIGNET)	0.2 mL	636	Dedham Neuro
SIG-39760-0500	ERAB (10D12) Monoclonal Antibody (SIGNET)	0.5 mL	1225	Dedham Neuro
SIG-39770-1000	ERAB (5F3) Monoclonal Antibody (SIGNET)	1.0 mL	2007	Dedham Neuro
SIG-39770-0200	ERAB (5F3) Monoclonal Antibody (SIGNET)	0.2 mL	636	Dedham Neuro
SIG-39770-0500	ERAB (5F3) Monoclonal Antibody (SIGNET)	0.5 mL	1225	Dedham Neuro
SIG-39780-1000	Apolipoprotein-D (GCDFP-24) Monoclonal Antibody (SIGNET)	1.0 mL	2007	Dedham Neuro
SIG-39780-0200	Apolipoprotein-D (GCDFP-24) Monoclonal Antibody (SIGNET)	0.2 mL	636	Dedham Neuro
SIG-39780-0500	Apolipoprotein-D (GCDFP-24) Monoclonal Antibody (SIGNET)	0.5 mL	1225	Dedham Neuro
SIG-39802-1000	Prion (4B4) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	2007	Dedham Neuro
SIG-39802-0200	Prion (4B4) Monoclonal Antibody, Purified (SIGNET)	0.2 mL	636	Dedham Neuro
SIG-39802-0500	Prion (4B4) Monoclonal Antibody, Purified (SIGNET)	0.5 mL	1225	Dedham Neuro
SIG-39805-1000	Prion (7D9) Monoclonal Antibody, Purified (SIGNET)	1.0 mL	2007	Dedham Neuro
SIG-39805-0200	Prion (7D9) Monoclonal Antibody, Purified (SIGNET)	0.2 mL	636	Dedham Neuro
SIG-39805-0500	Prion (7D9) Monoclonal Antibody, Purified (SIGNET)	0.5 mL	1225	Dedham Neuro
SIG-39810-0100	Prion (6D11) Monoclonal Antibody, Purified (SIGNET)	0.1 mL	581	Dedham Neuro
SIG-39810-0250	Prion (6D11) Monoclonal Antibody, Purified (SIGNET)	0.25 mL	1066	Dedham Neuro
SIG-39810-0500	Prion (6D11) Monoclonal Antibody, Purified (SIGNET)	0.5 mL	1828	Dedham Neuro
SIG-39830-1000	DJ-1 Monoclonal Antibody, Purified (SIGNET)	1.0 mL	1828	Dedham Neuro
SIG-39830-0200	DJ-1 Monoclonal Antibody, Purified (SIGNET)	0.2 mL	581	Dedham Neuro
SIG-39830-0500	DJ-1 Monoclonal Antibody, Purified (SIGNET)	0.5 mL	1115	Dedham Neuro
SIG-39835-050	DJ-1 Polyclonal Antibody, Purified (SIGNET)	0.05 mg	651	Dedham Neuro
SIG-39900-050	DJ-1 Human Protein Purified (SIGNET)	50 ug	582	Dedham Neuro
SIG-39938-020	sAPPbeta Recombinant Protein	0.02 mg	347	Dedham Neuro
SMI-21R-0100	GFAP (SMI 21) Monoclonal Antibody	0.1 mL	353	Covance Core

CatalogNo	Product Name	Size	EUR	Product Line
SMI-21R-0500	GFAP (SMI 21) Monoclonal Antibody	0.5 mL	534	Covance Core
SMI-22R-0100	GFAP (SMI 22) Monoclonal Antibody Cocktail	0.1 mL	353	Covance Core
SMI-22R-0500	GFAP (SMI 22) Monoclonal Antibody Cocktail	0.5 mL	534	Covance Core
SMI-23R-0100	GFAP (SMI 23) Monoclonal Antibody	0.1 mL	353	Covance Core
SMI-23R-0500	GFAP (SMI 23) Monoclonal Antibody	0.5 mL	534	Covance Core
SMI-24R-0100	GFAP (SMI 24) Monoclonal Antibody	0.1 mL	353	Covance Core
SMI-24R-0500	GFAP (SMI 24) Monoclonal Antibody	0.5 mL	534	Covance Core
SMI-25R-0100	GFAP (SMI 25) Monoclonal Antibody	0.1 mL	353	Covance Core
SMI-25R-0500	GFAP (SMI 25) Monoclonal Antibody	0.5 mL	534	Covance Core
SMI-26R-0100	GFAP (SMI 26) Monoclonal Antibody Cocktail	0.1 mL	353	Covance Core
SMI-26R-0500	GFAP (SMI 26) Monoclonal Antibody Cocktail	0.5 mL	534	Covance Core
SMI-310R-0100	Neurofilament H Phosphorylated (SMI 310) Monoclonal Antibody	0.1 mL	422	Covance Core
SMI-310R-0500	Neurofilament H Phosphorylated (SMI 310) Monoclonal Antibody	0.5 mL	733	Covance Core
SMI-311R-0100	Pan-Neuronal Neurofilament Marker (SMI 311) Monoclonal Antibody	0.1 mL	422	Covance Core
SMI-311R-0500	Pan-Neuronal Neurofilament Marker (SMI 311) Monoclonal Antibody	0.5 mL	733	Covance Core
SMI-312R-0100	Pan-Axonal Neurofilament Marker (SMI 312) Monoclonal Antibody	0.1 mL	422	Covance Core
SMI-312R-0500	Pan-Axonal Neurofilament Marker (SMI 312) Monoclonal Antibody	0.5 mL	733	Covance Core
SMI-31P-0100	Neurofilaments, Phosphorylated (SMI 31) Monoclonal Antibody, Purified	0,1 ml	476	Added April 10th
SMI-31R-0100	Neurofilaments, Phosphorylated (SMI 31) Monoclonal Antibody	0.1 mL	422	Covance Core
SMI-31R-0500	Neurofilaments, Phosphorylated (SMI 31) Monoclonal Antibody	0.5 mL	733	Covance Core
SMI-32P-0100	Neurofilament H Non-Phosphorylated (SMI 32) Monoclonal Antibody	0.1 mL	476	Covance Core
SMI-32R-0100	Neurofilament H Non-Phosphorylated (SMI 32) Monoclonal Antibody	0.1 mL	422	Covance Core
SMI-32R-0500	Neurofilament H Non-Phosphorylated (SMI 32) Monoclonal Antibody	0.5 mL	733	Covance Core
SMI-33R-0100	Neurofilament H & M Non-Phosphorylated (SMI 33) Monoclonal Antibody	0.1 mL	491	Covance Core
SMI-33R-0500	Neurofilament H & M Non-Phosphorylated (SMI 33) Monoclonal Antibody	0.5 mL	847	Covance Core
SMI-34R-0100	Neurofilament H Phosphorylated (SMI 34) Monoclonal Antibody	0.1 mL	422	Covance Core
SMI-34R-0500	Neurofilament H Phosphorylated (SMI 34) Monoclonal Antibody	0.5 mL	733	Covance Core
SMI-35R-0100	Neurofilaments, Hypophosphorylated (SMI 35) Monoclonal Antibody	0.1 mL	422	Covance Core
SMI-35R-0500	Neurofilaments, Hypophosphorylated (SMI 35) Monoclonal Antibody	0.5 mL	733	Covance Core
SMI-36R-0100	Neurofilament H Phosphorylated (SMI 36) Monoclonal Antibody	0.1 mL	422	Covance Core
SMI-36R-0500	Neurofilament H Phosphorylated (SMI 36) Monoclonal Antibody	0.5 mL	733	Covance Core
SMI-37R-0100	Neurofilament H Non-Phosphorylated (SMI 37) Monoclonal Antibody	0.1 mL	422	Covance Core
SMI-37R-0500	Neurofilament H Non-Phosphorylated (SMI 37) Monoclonal Antibody	0.5 mL	733	Covance Core
SMI-38R-0100	Neurofilament H Non-Phosphorylated (SMI 38) Monoclonal Antibody	0.1 mL	450	Covance Core
SMI-38R-0500	Neurofilament H Non-Phosphorylated (SMI 38) Monoclonal Antibody	0.5 mL	776	Covance Core
SMI-39R-0100	Neurofilament H Non-Phosphorylated (SMI 39) Monoclonal Antibody	0.1 mL	422	Covance Core
SMI-39R-0500	Neurofilament H Non-Phosphorylated (SMI 39) Monoclonal Antibody	0.5 mL	733	Covance Core
SMI-4010L-1000	ACTIVITY-SELECT Rabbit PAP	1.0 mL	160	Covance Core
SMI-4020L-1000	ACTIVITY-SELECT Goat PAP	1.0 mL	160	Covance Core
SMI-4050L-1000	Mouse ClonoPAP®	1.0 mL	298	Covance Core
SMI-4060L-1000	Rat PAP	1.0 mL	270	Covance Core
SMI-4070L-1000	ACTIVITY-SELECT Rat ClonoPAP®	1.0 mL	298	Covance Core
SMI-41R-0100	Gonadotropin-Releasing Hormone (LHRH) (SMI 41) Monoclonal Antibody	0.1 mL	353	Covance Core
SMI-41R-0500	Gonadotropin-Releasing Hormone (LHRH) (SMI 41) Monoclonal Antibody	0.5 mL	520	Covance Core
SMI-5010C-2000	Goat anti-mouse IgG	2.0 mL	69	Covance Core
SMI-5020C-2000	Rabbit anti-Goat IgG	2.0 mL	63	Covance Core
SMI-5030C-2000	Goat anti-Rabbit IgG	2.0 mL	63	Covance Core
SMI-5040C-2000	Goat anti-Rat IgG	2.0 mL	63	Covance Core
SMI-5050C-2000	Goat anti-Mouse IgG for Human Sections	2.0 mL	75	Covance Core
SMI-5060C-2000	Sheep anti-Rabbit IgG	2.0 mL	63	Covance Core
SMI-51R-0100	Tau (SMI 51) Monoclonal Antibody	0.1 mL	315	Covance Core
SMI-51R-0500	Tau (SMI 51) Monoclonal Antibody	0.5 mL	448	Covance Core
SMI-52R-0100	MAP2 (SMI 52) Monoclonal Antibody	0.1 mL	367	Covance Core
SMI-52R-0500	MAP2 (SMI 52) Monoclonal Antibody	0.5 mL	633	Covance Core
SMI-6010C-2000	Normal Rabbit Antibody	2.0 mL	56	Covance Core
SMI-6020C-2000	Normal Goat Antibody	2.0 mL	56	Covance Core
SMI-6040C-2000	Normal Rat Antibody	2.0 mL	63	Covance Core
SMI-6050C-2000	Normal Sheep Antibody	2.0 mL	56	Covance Core
SMI-61R-0500	Depolymerized β-Tubulin (SMI 61) Monoclonal Antibody	0.5 mL	392	Covance Core
SMI-62R-0500	Polymerized β-Tubulin (SMI 62) Monoclonal Antibody	0.5 mL	392	Covance Core
SMI-71R-0100	Rat Blood-Brain Barrier (SMI 71) Monoclonal Antibody	0.1 mL	353	Covance Core
SMI-71R-0500	Rat Blood-Brain Barrier (SMI 71) Monoclonal Antibody	0.5 mL	534	Covance Core
SMI-81R-0100	SNAP-25 (SMI 81) Monoclonal Antibody	0.1 mL	367	Covance Core
SMI-81R-0500	SNAP-25 (SMI 81) Monoclonal Antibody	0.5 mL	633	Covance Core
SMI-82R-0500	Synaptic Protein (SMI 82) Monoclonal Antibody	0.5 mL	362	Covance Core
SMI-91R-0100	Myelin CNPase (SMI 91) Monoclonal Antibody	0.1 mL	347	Covance Core
SMI-91R-0500	Myelin CNPase (SMI 91) Monoclonal Antibody	0.5 mL	591	Covance Core
SMI-94R-0100	Myelin Basic Protein (SMI 94) Monoclonal Antibody	0.1 mL	491	Covance Core
SMI-94R-0500	Myelin Basic Protein (SMI 94) Monoclonal Antibody	0.5 mL	847	Covance Core
SMI-99P-0100	Myelin Basic Protein (SMI 99) Monoclonal Antibody, Purified	0.1 mL	347	Covance Core
SMI-99P-0500	Myelin Basic Protein (SMI 99) Monoclonal Antibody, Purified	0.5 mL	591	Covance Core