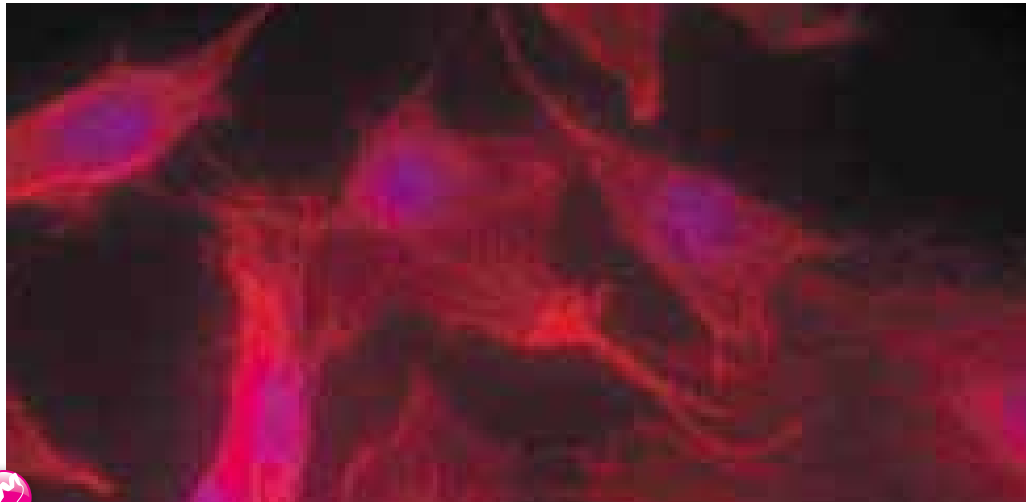


Assay Kits

The most extensive list of chromogenic and fluorogenic substrates for the sensitive analysis of enzyme activities. This validated and high quality set of products has been developed to perform your research on very diverse cellular processes.



The perfect combination between high quality peptides and the best dyes from the market

Benefits

- ▶ Ultra sensitive detection of enzyme activities
- ▶ ELISA of neurodegenerative disease-associated antibodies
- ▶ High-throughput screening of protease inhibitors in drug discovery

Peptides

Custom, Catalog & Labeled Peptides

Custom Peptides

- ▶ Long peptides (80-100 mer)
- ▶ Cyclic peptides
- ▶ Difficult sequences
- ▶ Large synthesis scale
- ▶ >95 % pure
- ▶ Stringent QC

Detection Reagents

HiLyte Fluor™ Dyes

- ▶ Proprietary fluorescent dyes
- ▶ Spanning the full visible spectrum
- ▶ More intense fluorescence than FITC, TAMRA and ROX
- ▶ More photostable and brighter than FITC, Cy® and DyLight®
- ▶ More soluble than other dyes

QXL™ Quenchers

- ▶ Proprietary quenchers
- ▶ Longer absorption wavelengths and higher extinction coefficients than DABCYL or DNP
- ▶ Individually optimized to pair with all popular fluorescent dyes

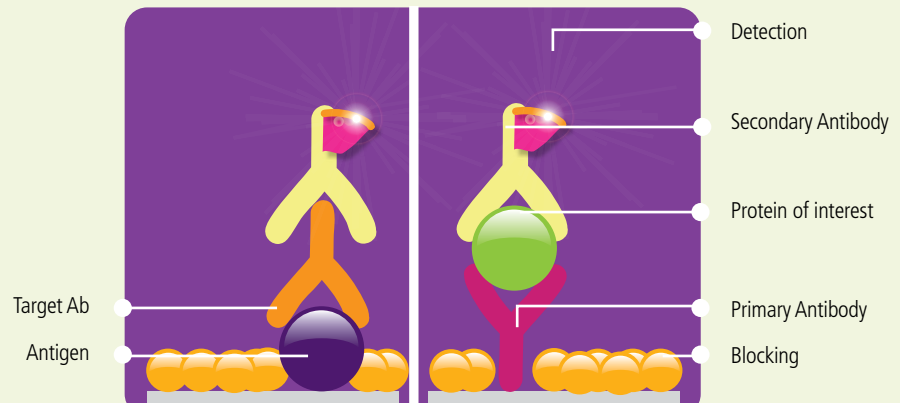
▶ Sensolyte® Assay Kits

- ▶ Large panel of Fluorescent / FRET peptides
- ▶ Custom Assay development

Detection methods used in Assay kits

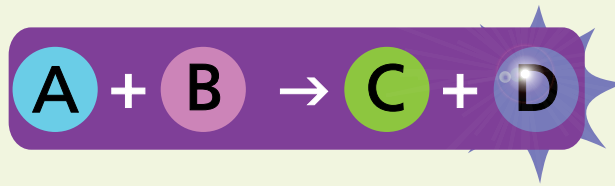
ELISA

During Enzyme-linked immunosorbent assay (ELISA), an unknown amount of target is specifically immobilized on a solid support (capture of target antibody by a specific antigen; capture of target protein by a specific antibody). Detection is performed using a labeled antibody conjugate, which produces a signal (fluorimetric, colorimetric or luminometric), proportional to the quantity of tested compound in the sample.



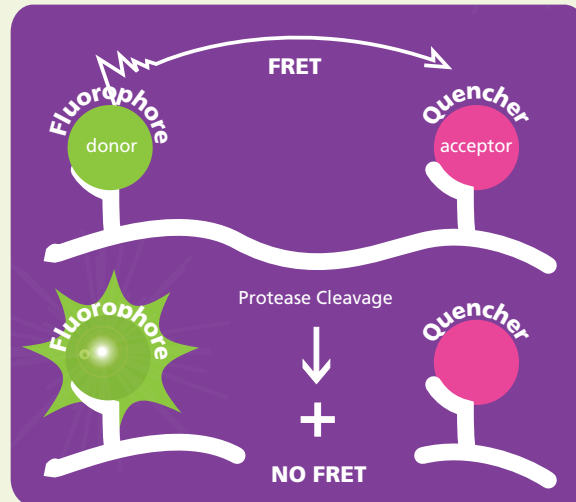
Enzymatic assay

All enzyme assays measure either the consumption of substrate or production of product over time. The concentration of substrates/products as well as the activity of enzymes can be assayed.



FRET protease assay

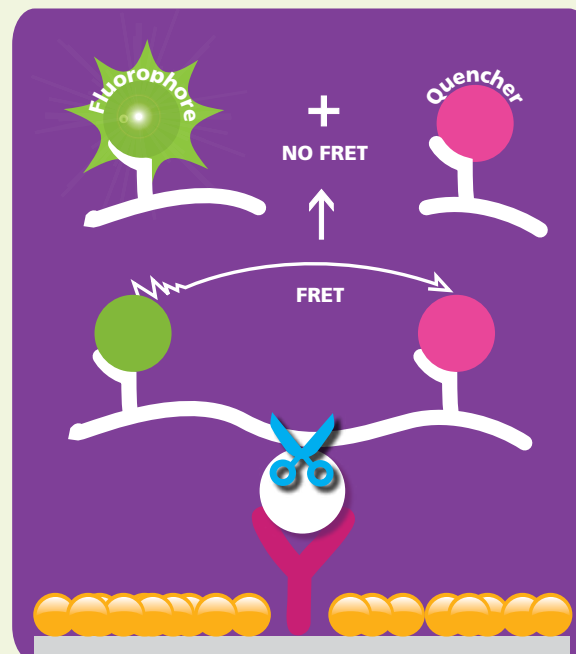
- Fluorescence or Förster resonance energy transfer (FRET) is a distance-dependent transfer of excited state energy from an excited donor (fluorophore) to an acceptor (another fluorophore or quencher). This transfer prevents the fluorophore to emit a fluorescent signal at its emission wavelength. Consequently these assays measure events allowing the physical separation of the fluorophore and the quencher, such as proteolytic cleavage, leading to the release of the fluorescent signal.
- Fluorogenic peptide protease assays also measure the activity of proteases but a fluorogenic peptide is used as a substrate.



FRET + ELISA

(Sensolyte® Plus kits)

FRET technology combined with ELISA assay allows a specific detection of the activity of a particular protease in a biological sample, possibly containing multiple proteases which recognize the same FRET peptide-substrate. The target protease is selectively captured by ELISA, then its activity is assayed using a FRET peptide.



Assay kit overview

Enzyme/Antibody	Product Detection	Reference
Angiotensin-converting enzyme 2 (ACE2)	SensoLyte® 390 ACE2 Activity Assay Kit - 100 assays - Fluorimetric	AS-72086
Aggrecanase-1	SensoLyte® 520 Aggrecanase-1 Assay Kit - 100 assays - Fluorimetric	AS-72114
α-Secretase	SensoLyte® 520 TACE (α-Secretase) Activity Assay Kit - 100 assays - Fluorimetric	AS-72085
α-Synuclein	SensoLyte® Anti - α - Synuclein (Human, Mouse, Rat) ELISA Kit	AS-55550
β-Galactosidase	SensoLyte® FDG β-Galactosidase Assay Kit - 500 assays - Fluorimetric	AS-72133
	SensoLyte® MUG β-Galactosidase Assay Kit - 500 assays - Fluorimetric	AS-72132
	SensoLyte® ONPG β-Galactosidase Assay Kit - 500 assays - Colorimetric	AS-72134
Biotin	SensoLyte® Biotin Quantitation Kit - 200 assays - Fluorimetric	AS-72163
	SensoLyte® Calcein Cell Viability Assay Kit - 1000 assays - Fluorimetric	AS-72126
Cell Proliferation & Cytotoxicity	DHL™ Cell Viability and Proliferation Assay Kit - 2000 assays - Fluorimetric	AS-71300
	DHL™ Cell Viability and Proliferation Assay Kit - 10000 assays - Fluorimetric	AS-71301
	DHL™ Cell Cytotoxicity Assay Kit - 500 assays - Fluorimetric	AS-71302
	DHL™ Cell Cytotoxicity Assay Kit - 5000 assays - Fluorimetric	AS-71303
Calpain	SensoLyte® 520 Calpain Activity Assay Kit - 100 assays - Fluorimetric	AS-72149
	SensoLyte® AMC Calpain Activity Assay Kit - 100 assays - Fluorimetric	AS-72150
Caspases	SensoLyte® AFC Caspase Profiling Kit - 2 x 96-well plate - Fluorimetric	AS-71116
	SensoLyte® AMC Caspase Profiling Kit - 2 x 96-well plate - Fluorimetric	AS-71120
	SensoLyte® AFC Caspase Substrate Sampler Kit - 8 x 100 assays - Fluorimetric	AS-71117
	SensoLyte® AMC Caspase Substrate Sampler Kit - 8 x 100 assays - Fluorimetric	AS-71121
	SensoLyte® Homogeneous AnaRed Caspase-3/7 Assay Kit - 500 assays - Fluorimetric	AS-71122
	SensoLyte® Homogeneous Rh110 Caspase-3/7 Assay Kit - 500 assays - Fluorimetric	AS-71141
	SensoLyte® Homogeneous AFC Caspase-3/7 Assay Kit - 500 assays - Fluorimetric	AS-71114
	SensoLyte® Homogeneous AMC Caspase-3/7 Assay Kit - 500 assays - Fluorimetric	AS-71118
	SensoLyte® Homogeneous AFC Caspase - 8 Assay Kit - 100 assays - Fluorimetric	AS-72088-100
	SensoLyte® Homogeneous AFC Caspase - 8 Assay Kit - 200 assays - Fluorimetric	AS-72088-200
Cathepsins	SensoLyte® 520 Cathepsin B Assay Kit - 100 assays - Fluorimetric	AS-72164
	SensoLyte® 440 Cathepsin B Assay Kit - 100 assays - Fluorimetric	AS-72165
	SensoLyte® 520 Cathepsin D Assay Kit - 100 assays - Fluorimetric	AS-72170
	SensoLyte® 390 Cathepsin D Assay Kit - 100 assays - Fluorimetric	AS-72098
	SensoLyte® 520 Cathepsin G Assay Kit - 100 assays - Fluorimetric	AS-72185
	SensoLyte® Rh110 Cathepsin G Assay Kit - 100 assays - Fluorimetric	AS-72186
	SensoLyte® Rh110 Cathepsin K Assay Kit - 100 assays - Fluorimetric	AS-72152
	SensoLyte® 520 Cathepsin K Assay Kit - 100 assays - Fluorimetric	AS-72171
	SensoLyte® 520 Cathepsin S Assay Kit - 100 assays - Fluorimetric	AS-72099
	SensoLyte® 440 Cathepsin S Assay Kit - 100 assays - Fluorimetric	AS-72100
Deubiquitination	SensoLyte® 520 Deubiquitination Assay Kit - 100 assays - Fluorimetric	AS-72203
	SensoLyte® 440 Deubiquitination Assay Kit - 100 assays - Fluorimetric	AS-72204
DPPIV	SensoLyte® Rh110 DPPIV Assay Kit - 100 assays - Fluorimetric	AS-72196
	SensoLyte® AMC DPPIV Assay Kit - 100 assays - Fluorimetric	AS-72197
Elastase	SensoLyte® Green Elastase Assay Kit - 500 assays - Fluorimetric	AS-72178
	SensoLyte® Rh110 Elastase Assay Kit - 100 assays - Fluorimetric	AS-72179
Factor Xa	SensoLyte® Rh110 Factor Xa Assay Kit - Fluorimetric	AS-72207
	SensoLyte® 520 Factor Xa Assay Kit - Fluorimetric	AS-72208
Generic Protease	SensoLyte® Green Protease Assay Kit - 500 assays - Fluorimetric	AS-71124
	SensoLyte® Red Protease Assay Kit - 500 assays - Fluorimetric	AS-71140
Glutathione/ Glutathione -S-Transferase	SensoLyte® Total GSH Assay Kit - 500 assays - Colorimetric	AS-72153
	SensoLyte® 520 Total GSH Assay Kit - 100 assays - Fluorimetric	AS-72154
	SensoLyte® GST Activity Assay Kit - 100 assays - Fluorimetric	AS-72157
HCV	SensoLyte® Glutathione Cellular Assay Kit - 100 assays - Fluorimetric	AS-72158
	SensoLyte® 620 HCV Protease Assay Kit - 100 assays - Fluorimetric	AS-71146
	SensoLyte® 520 HCV Protease Assay Kit - 100 assays - Fluorimetric	AS-71145
	SensoLyte® 490 HCV Protease Assay Kit - 500 assays - Fluorimetric	AS-71126
	SensoLyte® 490 HCV Protease Assay Kit - 200 assays - Fluorimetric	AS-72087
Histone Acetyl Transferases (HAT)	SensoLyte® HAT (p300) Assay Kit - 100 assays - Fluorimetric	AS-72172
	SensoLyte® HAT (pCAF) Assay Kit - 100 assays - Fluorimetric	AS-72173
Histone Sampler kits	SensoLyte® Histone H3 (1-21) Sampler Kit - Biotinylated	AS-72175
	SensoLyte® Methylated Histone H3 (1-21) Sampler Kit - Biotinylated	AS-72176
Histone Deacetylases (HDAC)	SensoLyte® 520 HDAC Activity Assay Kit - 100 assays - Fluorimetric	AS-72084
	SensoLyte® 440 HDAC Activity Assay Kit - 100 assays - Fluorimetric	AS-72083
Sirtuin1 (Sirt1)	SensoLyte® 520 FRET Sirt1 Assay Kit - 100 assays - Fluorimetric	AS-72155
	SensoLyte® Green Sirt1 Assay Kit - 100 assays - Fluorimetric	AS-72156
Sirtuin2 (Sirt2)	SensoLyte® Green Sirt2 Assay Kit - 100 assays - Fluorimetric	AS-72188
	SensoLyte® 520 FRET Sirt2 Assay Kit - 100 assays - Fluorimetric	AS-72189
Histone Demethylases (LSD1)	SensoLyte® LSD1 Assay Kit - 100 assays - Fluorimetric	AS-72131
HIV	SensoLyte® 620 HIV Protease Assay Kit - 100 assays - Fluorimetric	AS-71148
	SensoLyte® 520 HIV Protease Assay Kit - 100 assays - Fluorimetric	AS-71147
	SensoLyte® 490 HIV Protease Assay Kit - 500 assays - Fluorimetric	AS-71127
Kinases	SensoLyte® Protein Kinase Substrate Sampler Kit - Biotinylated	AS-71200
	SensoLyte® Serine/Threonine Protein Kinase Substrate Sampler Kit -1 - FAM-labeled	AS-71201
	SensoLyte® Serine/Threonine Protein Kinase Substrate Sampler Kit-2 - FAM-Labeled	AS-71202
	SensoLyte® Serine/Threonine Protein Kinase Substrate Sampler Kit-1 - TAMRA-labeled	AS-72029
	SensoLyte® Serine/Threonine Protein Kinase Substrate Sampler Kit-2 -TAMRA-labeled	AS-72030
	SensoLyte® Tyrosine Protein Kinase Substrate Profiling Kit - Fluorimetric	AS-71203
Matrix Metalloproteinases (MMPs)	SensoLyte® 570 Generic MMP Assay Kit - 100 assays - Fluorimetric	AS-72101
	SensoLyte® 520 Generic MMP Activity Kit - 100 assays - Fluorimetric	AS-71158
	SensoLyte® Generic MMP Assay Kit - 100 assays - Colorimetric	AS-72095
	SensoLyte® 390 Generic MMP Activity Kit - 100 assays - Fluorimetric	AS-72202
	SensoLyte® 520 MMP Profiling Kit -2 x 96-well plate - Fluorimetric	AS-71136
MMP-1	SensoLyte® 520 MMP Substrate Sampler Kit - 16 x 20 assays - Fluorimetric	AS-71170
	SensoLyte® Plus 520 MMP-1 Assay Kit - 96-well plate - Fluorimetric and Enhanced Selectivity	AS-72012
	SensoLyte® 520 MMP-1 Assay Kit - 100 assays - Fluorimetric	AS-71150
	SensoLyte® 490 MMP-1 Assay Kit - 500 assays - Fluorimetric	AS-71128
MMP-2	SensoLyte® MMP-1 ELISA Kit - 96-well plate - Colorimetric	AS-72102
	SensoLyte® 520 MMP-2 Assay Kit - 100 assays - Fluorimetric	AS-71151
	SensoLyte® 490 MMP-2 Assay Kit - 500 assays - Fluorimetric	AS-71129
	SensoLyte® MMP-2 ELISA Kit - 96-well plate - Colorimetric	AS-72148

Enzyme/Antibody	Product Detection	Reference
MMP-3	Sensolyte® 520 MMP-3 Assay Kit - 100 assays - Fluorimetric	AS-71152
	Sensolyte® 490 MMP-3 Assay Kit - 500 assays - Fluorimetric	AS-71130
	Sensolyte® MMP-3 ELISA Kit - 96-well plate - Colorimetric	AS-72103
MMP-7	Sensolyte® 520 MMP-7 Assay Kit - 100 assays - Fluorimetric	AS-71153
	Sensolyte® 490 MMP-7 Assay Kit - 500 assays - Fluorimetric	AS-71132
MMP-8	Sensolyte® 520 MMP-8 Assay Kit - 100 assays - Fluorimetric	AS-71154
	Sensolyte® 490 MMP-8 Assay Kit - 500 assays - Fluorimetric	AS-71133
	Sensolyte® MMP-8 ELISA Kit - 96-well plate - Colorimetric	AS-72104
MMP-9	Sensolyte® Plus 520 MMP-9 Assay Kit - 96-well plate - Fluorimetric and Enhanced Selectivity	AS-72017
	Sensolyte® 520 MMP-9 Assay Kit - 100 assays - Fluorimetric	AS-71155
	Sensolyte® 490 MMP-9 Assay Kit - 500 assays - Fluorimetric	AS-71134
	Sensolyte® MMP-9 ELISA Kit - 96-well plate - Colorimetric	AS-72105
MMP-10	Sensolyte® 520 MMP-10 Assay Kit - 100 assays - Fluorimetric	AS-72024
	Sensolyte® MMP-10 ELISA Kit - 96-well plate - Colorimetric	AS-72106
MMP-12	Sensolyte® 520 MMP-12 Assay Kit - 100 assays - Fluorimetric	AS-71157
	Sensolyte® 490 MMP-12 Assay Kit - 500 assays - Fluorimetric	AS-71137
MMP-13	Sensolyte® Plus 520 MMP-13 Assay Kit - 96-well plate - Fluorimetric and Enhanced Selectivity	AS-72019
	Sensolyte® 520 MMP-13 Assay Kit - 100 assays - Fluorimetric	AS-71156
	Sensolyte® 490 MMP-13 Assay Kit - 500 assays - Fluorimetric	AS-71135
	Sensolyte® MMP-13 ELISA Kit - 96-well plate - Colorimetric	AS-72107
MMP-14	Sensolyte® 520 MMP-14 Assay Kit - 100 assays - Fluorimetric	AS-72025
MMP-15	Sensolyte® 520 MMP-15 Assay Kit - 100 assays - Fluorimetric	AS-72198
MMP-16	Sensolyte® 520 MMP-16 Assay Kit - 100 assays - Fluorimetric	AS-72200
MMP-24	Sensolyte® 520 MMP-24 Assay Kit - 100 assays - Fluorimetric	AS-72201
Myelin Oligodendrocyte Glycoprotein (MOG)	Sensolyte® Anti-Human MOG (1-125) IgG Quantitative ELISA Kit - 12 x 8 well strips - Colorimetric	AS-55153
	Sensolyte® Anti-Mouse MOG (1-125) IgG Quantitative ELISA Kit - 12 x 8 well strips - Colorimetric	AS-55156
	Sensolyte® Anti-Rat MOG (1-125) IgG Quantitative ELISA Kit - 12 x 8 well strips - Colorimetric	AS-55157
	Sensolyte® Anti-MOG (35-55) IgG Quantitative ELISA Kit (mouse/rat) - 12 x 8 well strips - Colorimetric	AS-54465
NADP/NADPH	Sensolyte® NADP/NADPH Assay Kit - 100 assays - Colorimetric	AS-72205
	Sensolyte® NADP/NADPH Assay Kit - 100 assays - Fluorimetric	AS-72206
Peroxidase/ Hydrogen Peroxyde	Sensolyte® ADHP Peroxidase ELISA Assay Kit (mouse) - 500 assays - Fluorimetric	AS-71110-M
	Sensolyte® ADHP Peroxidase ELISA Assay Kit (rat) - 500 assays - Fluorimetric	AS-71110-R
	Sensolyte® Luminescent Peroxidase ELISA Assay Kit - 500 assays - Luminometric	AS-72128
	Sensolyte® Luminescent Peroxidase Assay Kit - 500 assays - Luminometric	AS-72127
	Sensolyte® ADHP Hydrogen Peroxide Assay Kit - 500 assays - Fluorimetric	AS-71112
Phosphatase	Sensolyte® MG Phosphate Assay Kit - 1000 assays - Colorimetric	AS-71103
	Sensolyte® pNPP Protein Phosphatase Assay Kit - 500 assays - Colorimetric	AS-71105
	Sensolyte® FDP Protein Phosphatase Assay Kit - 500 assays - Fluorimetric	AS-71100
	Sensolyte® MFP Protein Phosphatase Assay Kit - 500 assays - Fluorimetric	AS-71104
	Sensolyte® MFP Acid Phosphatase Assay Kit - 500 assays - Fluorimetric	AS-71231
	Sensolyte® pNPP Alkaline Phosphatase Assay Kit - 500 assays - Colorimetric	AS-72146
Alkaline Phosphatase	Sensolyte® FDP Alkaline Phosphatase Assay Kit - 500 assays - Fluorimetric	AS-71109
	Sensolyte® Luminescent Alkaline Phosphatase Assay Kit - 500 assays - Luminometric	AS-72122
	Sensolyte® FDP Alkaline Phosphatase ELISA Assay Kit - 500 assays - Fluorimetric	AS-71101-M
	Sensolyte® FDP Alkaline Phosphatase ELISA Assay Kit - 500 assays - Fluorimetric	AS-71101-R
	Sensolyte® pNPP Alkaline Phosphatase ELISA Assay Kit - 500 assays - Colorimetric	AS-72147-G
	Sensolyte® pNPP Alkaline Phosphatase ELISA Assay Kit - 500 assays - Colorimetric	AS-72147-M
	Sensolyte® pNPP Alkaline Phosphatase ELISA Assay Kit - 500 assays - Colorimetric	AS-72147-R
	Sensolyte® Luminescent Alkaline Phosphatase ELISA Assay Kit - 500 assays - Luminometric	AS-72123
	Sensolyte® pNPP Secreted Alkaline Phosphatase Reporter Gene Assay Kit - 500 assays - Colorimetric	AS-72144
	Sensolyte® FDP Secreted Alkaline Phosphatase Reporter Gene Assay Kit - 500 assays - Fluorimetric	AS-71108
Sensolyte® Luminescent Secreted Alkaline Phosphatase Reporter Gene Assay Kit - 500 assays - Luminometric	AS-72145	
Plasmin	Sensolyte® Rh110 Plasmin Activity Assay Kit - 100 assays - Fluorimetric	AS-72125
	Sensolyte® AFC Plasmin Activity Assay Kit - 100 assays - Fluorimetric	AS-72124
Plasminogen Activator	Sensolyte® AFC Urokinase (uPA) Activity Assay Kit - 100 assays - Fluorimetric	AS-72159
Protein Quantitation (OPA)	Sensolyte® AMC-Tissue Type (tPA) Activity Assay Kit - 100 assays - Fluorimetric	AS-72160
Proteolipid Protein (PLP)	Sensolyte® OPA Protein Quantitation Kit - 500 assays - Fluorimetric	AS-71015
	Sensolyte® Anti-PLP (139-151) IgG Quantitative ELISA Kit (mouse) - 12 x 8 well strips	AS-55524
Renin	Sensolyte® Anti-PLP (178-191) IgG Quantitative ELISA Kit (mouse) - 12 x 8 well strips	AS-55523
	Sensolyte® 520 Rat Renin Assay Kit - 100 assays - Fluorimetric	AS-72140
	Sensolyte® 520 Renin Assay Kit - 100 assays - Fluorimetric	AS-72040
	Sensolyte® 390 Renin Assay Kit - 100 assays - Fluorimetric	AS-72039
	Sensolyte® 520 Mouse Renin Assay Kit - 100 assays - Fluorimetric	AS-72161
Thiol	Sensolyte® Thiol Quantitation Assay Kit - 500 assays - Colorimetric	AS-72136
	Sensolyte® ABD-F Thiol Quantitation Assay Kit - 500 assays - Fluorimetric	AS-72137
	Sensolyte® 520 Thiol Quantitation Assay Kit - 500 assays - Fluorimetric	AS-72138
Thrombin	Sensolyte® 520 Thrombin Activity Assay Kit - Fluorimetric	AS-72129
	Sensolyte® AFC Thrombin Activity Assay Kit - Fluorimetric	AS-72130
WNV (West Nile Virus)	Sensolyte® 570 West Nile Virus Protease Assay Kit - Fluorimetric	AS-72080
	Sensolyte® 440 West Nile Virus Protease Assay Kit - Fluorimetric	AS-72079

Cy® is a registered trademark of GE Healthcare Bio-Sciences Corp.

DHL™ is a trademark of AnaSpec, Inc.

Dylight® is a registered trademark of Thermo Fischer Scientific Inc and its subsidiaries.

HiLyte Fluor™ is a trademark of AnaSpec, Inc.

QXL™ is a trademark of AnaSpec, Inc.

Sensolyte® is a registered trademark of AnaSpec, Inc.

License Statement

Cy⁵ amidite, Cy³ amidite, Cy^{3.5} amidite, and Cy^{5.5} amidite

The purchase of products with or in which Cy⁵ amidite, Cy³ amidite, Cy^{3.5} amidite, and Cy^{5.5} amidite may be used or combined, includes a limited non-exclusive sublicense under U.S. Patents 5,556,959 and 5,808,044 and foreign equivalent patents and other foreign and U.S. counterpart applications to use the amidites in such product to perform research. NO OTHER LICENSE IS GRANTED EXPRESSLY, IMPLIEDLY OR BY ESTOPPEL. Use of such product for commercial purposes is strictly prohibited without written permission from Amersham Biosciences Corp. For information concerning availability of additional licenses to

practise the patented methodologies, please contact Amersham Biosciences Corp., Business Licensing Manager, Amersham Place, Little Chalfont, Bucks, HP79NA.

HiLyte™ dyes

Use of this product employs the following patent rights licensed to the purchaser from AnaSpec, Inc.: (a) the claims of 7,465,810; 7,754,893; 7,820,783; 7,790,394; 7,985,552; 7,989,638 (b) any claims issuing from U.S. Patent Applications Serial No. 12/287,074; 12/583,513; 12/925,505 and (c) all patents to be issued pursuant thereto, and all continuations, continuations-in-part, reissues, substitutes, and extensions thereof. The

purchase of this product includes a limited, non-transferable immunity from suit under the foregoing patent claims for using only this amount of product for the purchaser's own internal research. No right under any other patent claim, no right to perform any patented method, and no right to perform commercial services of any kind, including without limitation reporting the results of purchaser's activities for a fee or other commercial consideration, is conveyed expressly, by implication, or by estoppel. This product is for research use only. Diagnostic uses under AnaSpec, Inc. patents require a separate license from AnaSpec, Inc.

More information

How to order?

www.eurogentec.com

order@eurogentec.com

+32 4 264 07 88

Need support?

info@eurogentec.com

00 800 666 00 123

[European Toll Free Number]

Eurogentec
Experiences true partnership

ANA SPEC