

# SensoLyte<sup>®</sup> Assay Kits

Detailed protocols • Convenient 96-well formats Ideal for drug discovery & HTS

Trusted by researchers worldwide for over **two decades**, and featured in high impact journals, **SensoLyte<sup>®</sup> Assay Kits** offer **accurate, sensitive, and user-friendly** detection, quantification and/or activity measurement of various **enzymes, analytes, reporter genes** and more. Most kits utilize **fluorescent, colorimetric, or luminescent** detection methods.

## Enzyme Activity Assay kits

Target	Catalog #	Product name	Detection	Color
3CL Protease	AS-72262	SensoLyte <sup>®</sup> 520 SARS-CoV-2 3CL Protease Activity Assay Kit	Fluorimetric	Green
Acetylcholinesterase	AS-72242	SensoLyte <sup>®</sup> 520 Acetylcholinesterase Activity Assay Kit	Fluorimetric	Green
ADAM10	AS-72226	SensoLyte <sup>®</sup> 520 ADAM10 Activity Assay Kit	Fluorimetric	Green
ADAMTS-13	AS-72232	SensoLyte <sup>®</sup> 520 ADAMTS13 Activity Assay Kit	Fluorimetric	Green
ADAMTS-4 (Aggrecanase)	AS-72114	SensoLyte <sup>®</sup> 520 Aggrecanase-1 Activity Assay Kit	Fluorimetric	Green
Secretase	AS-72085	SensoLyte <sup>®</sup> 520 TACE ( $\alpha$ -Secretase) Activity Assay Kit	Fluorimetric	Green
	AS-71144	SensoLyte <sup>®</sup> 520 b-Secretase (BACE1) Activity Assay Kit	Fluorimetric	Green
	AS-72225	SensoLyte <sup>®</sup> 520 BACE2 Activity Assay Kit	Fluorimetric	Green
Angiotensin Converting Enzyme	AS-72086	SensoLyte <sup>®</sup> 390 ACE2 Activity Assay Kit	Fluorimetric	Blue
Calpain	AS-72149	SensoLyte <sup>®</sup> 520 Calpain Activity Assay Kit	Fluorimetric	Green
	AS-72150	SensoLyte <sup>®</sup> AMC Calpain Activity Assay Kit	Fluorimetric	Blue
Caspase	AS-71114	SensoLyte <sup>®</sup> Homogeneous AFC Caspase-3/7 Activity Assay Kit	Fluorimetric	Blue
	AS-71118	SensoLyte <sup>®</sup> Homogeneous AMC Caspase-3/7 Activity Assay Kit	Fluorimetric	Blue
	AS-71141	SensoLyte <sup>®</sup> Homogeneous Rh110 Caspase-3/7 Activity Assay Kit	Fluorimetric	Green
Cathepsin	AS-72165	SensoLyte <sup>®</sup> 440 Cathepsin B Activity Assay Kit	Fluorimetric	Blue
	AS-72164	SensoLyte <sup>®</sup> 520 Cathepsin B Activity Assay Kit	Fluorimetric	Green
	AS-72098	SensoLyte <sup>®</sup> 390 Cathepsin D Activity Assay Kit	Fluorimetric	Blue
	AS-72170	SensoLyte <sup>®</sup> 520 Cathepsin D Activity Assay Kit	Fluorimetric	Green
	AS-72222	SensoLyte <sup>®</sup> 520 Cathepsin E Activity Assay Kit	Fluorimetric	Green
	AS-72185	SensoLyte <sup>®</sup> 520 Cathepsin G Activity Assay Kit	Fluorimetric	Green
	AS-72171	SensoLyte <sup>®</sup> 520 Cathepsin K Activity Assay Kit	Fluorimetric	Green
	AS-72218-	SensoLyte <sup>®</sup> 520 Cathepsin L Activity Assay Kit	Fluorimetric	Green
AS-72099	SensoLyte <sup>®</sup> 520 Cathepsin S Activity Assay Kit	Fluorimetric	Green	
Deubiquitinating enzyme	AS-72203	SensoLyte <sup>®</sup> 520 Deubiquitination Assay Kit	Fluorimetric	Green
Dipeptidylaminopeptidase (DPPIV)	AS-72196-	SensoLyte <sup>®</sup> Rh110 DPPIV Activity Assay Kit	Fluorimetric	Green
	AS-72197-	SensoLyte <sup>®</sup> AMC DPPIV Activity Assay Kit	Fluorimetric	Blue

# Enzyme Activity Assay kits

Target	Catalog #	Product name	Detection	Color
Elastase	AS-72178	SensoLyte® Green Elastase Activity Assay Kit	Fluorimetric	Green
	AS-72179	SensoLyte® Rh110 Elastase Activity Assay Kit	Fluorimetric	Green
Endothelin Converting Enzyme (ECE)	AS-72243	SensoLyte® 520 ECEs Activity Assay Kit	Fluorimetric	Green
Enterokinase	AS-72209	SensoLyte® 520 Enterokinase Activity Assay Kit	Fluorimetric	Green
	AS-72210	SensoLyte® Rh110 Enterokinase Activity Assay Kit	Fluorimetric	Green
Factor Xa	AS-72208	SensoLyte® 520 Factor Xa Activity Assay Kit	Fluorimetric	Green
	AS-72207	SensoLyte® Rh110 Factor Xa Activity Assay Kit	Fluorimetric	Green
Furin	AS-72256	SensoLyte® Rh110 Furin Activity Assay Kit	Fluorimetric	Green
Glucocerebrosidase	AS-72258	SensoLyte® Blue Glucocerebrosidase (GBA) Activity Assay Kit	Fluorimetric	Blue
	AS-72259	SensoLyte® Red Glucocerebrosidase (GBA) Activity Assay Kit	Fluorimetric	Orange
Glutaminyl cyclase	AS-72230	SensoLyte® Green Glutaminyl Cyclase Activity Assay Kit	Fluorimetric	Green
Granzyme A	AS-72260	SensoLyte® 520 Granzyme A Activity Assay Kit	Fluorimetric	Green
Granzyme B	AS-72261	SensoLyte® 520 Granzyme B Activity Assay Kit	Fluorimetric	Green
Glutathione S Transferase (GST)	AS-72157	SensoLyte® GST Activity Assay Kit	Fluorimetric	Blue
Hepatitis C Virus (HCV) Protease	AS-71126	SensoLyte® 490 HCV Protease Activity Assay Kit - 500 assays	Fluorimetric	Blue
	AS-71145	SensoLyte® 520 HCV Protease Activity Assay Kit	Fluorimetric	Green
Histone acetyltransferase (HAT)	AS-72172	SensoLyte® HAT (p300) Activity Assay Kit	Fluorimetric	Blue
Histone deacetylase (HDAC)	AS-72084	SensoLyte® 520 HDAC Activity Assay Kit	Fluorimetric	Green
Human Immunodeficiency Virus 1 (HIV1) protease	AS-71127	SensoLyte® 490 HIV Protease Activity Assay Kit	Fluorimetric	Blue
	AS-71147	SensoLyte® 520 HIV Protease Activity Assay Kit	Fluorimetric	Green
Insulin Degrading Enzyme (IDE)	AS-72231	SensoLyte® 520 IDE Activity Assay Kit	Fluorimetric	Green
Kallikrein	AS-72255	SensoLyte® Rh110 Plasma Kallikrein Activity Assay Kit	Fluorimetric	Green
Matriptase	AS-72241	SensoLyte® Rh110 Matriptase Activity Assay Kit	Fluorimetric	Green
Matrix Metalloproteinase (MMP) Generic	AS-71158	SensoLyte® 520 Generic MMP Activity Assay Kit	Fluorimetric	Green
	AS-71136	SensoLyte® 520 MMP Profiling Kit	Fluorimetric	Green
	AS-71170	SensoLyte® 520 MMP Substrate Sampler Kit	Fluorimetric	Green
	AS-72095	SensoLyte® Generic MMP Activity Assay Kit	Colorimetric	n/a
	AS-72101	SensoLyte® 570 Generic MMP Activity Assay Kit	Fluorimetric	Orange
MMP 1, Interstitial collagenase	AS-71128	SensoLyte® 490 MMP-1 Activity Assay Kit	Fluorimetric	Blue
	AS-71150	SensoLyte® 520 MMP-1 Activity Assay Kit	Fluorimetric	Green
	AS-72012	SensoLyte® Plus 520 MMP-1 Activity Assay Kit - Enhanced Selectivity	Fluorimetric	Green
MMP 2, Gelatinase-A	AS-71129	SensoLyte® 490 MMP-2 Activity Assay Kit	Fluorimetric	Blue
	AS-71151	SensoLyte® 520 MMP-2 Activity Assay Kit	Fluorimetric	Green
MMP 3, Stromelysin 1	AS-71130	SensoLyte® 490 MMP-3 Activity Assay Kit	Fluorimetric	Blue
	AS-71152	SensoLyte® 520 MMP-3 Activity Assay Kit	Fluorimetric	Green
MMP 7, Matrilysin	AS-71153	SensoLyte® 520 MMP-7 Activity Assay Kit	Fluorimetric	Green
MMP 8, Neutrophil collagenase	AS-71133	SensoLyte® 490 MMP-8 Activity Assay Kit	Fluorimetric	Blue
	AS-71154	SensoLyte® 520 MMP-8 Activity Assay Kit	Fluorimetric	Green
MMP 9, Gelatinase-B	AS-71134	SensoLyte® 490 MMP-9 Activity Assay Kit	Fluorimetric	Blue
	AS-71155	SensoLyte® 520 MMP-9 Activity Assay Kit	Fluorimetric	Green
	AS-72017	SensoLyte® Plus 520 MMP-9 Activity Assay Kit - Enhanced Selectivity	Fluorimetric	Green
MMP 10, Stromelysin 2	AS-72024	SensoLyte® 520 MMP-10 Activity Assay Kit	Fluorimetric	Green
MMP 12, Macrophage metalloelastase	AS-71137	SensoLyte® 490 MMP-12 Activity Assay Kit	Fluorimetric	Blue
	AS-71157	SensoLyte® 520 MMP-12 Activity Assay Kit	Fluorimetric	Green
MMP 13, Collagenase 3	AS-71135	SensoLyte® 490 MMP-13 Activity Assay Kit	Fluorimetric	Blue
	AS-71156	SensoLyte® 520 MMP-13 Activity Assay Kit	Fluorimetric	Green
	AS-72019	SensoLyte® Plus 520 MMP-13 Activity Assay Kit - Enhanced Selectivity	Fluorimetric	Green
MMP 14, MT1-MMP	AS-72025	SensoLyte® 520 MMP-14 Activity Assay Kit	Fluorimetric	Green

Target	Catalog #	Product name	Detection	Color
Meprin	AS-72253	SensoLyte® 520 Meprin $\alpha$ Activity Assay Kit	Fluorimetric	Green
	AS-72254	SensoLyte® 520 Meprin $\beta$ Activity Assay Kit	Fluorimetric	Green
Nepriylsin	AS-72223	SensoLyte® 520 Nepriylsin Activity Assay Kit	Fluorimetric	Green
Papain-like Protease (SARS CoV 2 PLPro)	AS-72274	SensoLyte® 520 SARS-CoV-2 Papain-like Protease/Deubiquitinase Activity Assay Kit	Fluorimetric	Green
Peptidyl-prolyl isomerase (Pini)	AS-72240	SensoLyte® Green PINI Activity Assay Kit	Fluorimetric	Green
Plasmin	AS-72124	SensoLyte® AFC Plasmin Activity Assay Kit	Fluorimetric	Blue
	AS-72125	SensoLyte® Rh110 Plasmin Activity Assay Kit	Fluorimetric	Green
Plasminogen activator	AS-72159	SensoLyte® AFC Urokinase (uPA) Activity Assay Kit	Fluorimetric	Blue
	AS-72160	SensoLyte® AMC Tissue-type (tPA) Activity Assay Kit	Fluorimetric	Blue
Phosphatase	AS-71100	SensoLyte® FDP Protein Phosphatase Assay Kit	Fluorimetric	Green
	AS-71105	SensoLyte® pNPP Protein Phosphatase Assay Kit	Colorimetric	n/a
Protease (generic)	AS-71124	SensoLyte® Green Protease Assay Kit	Fluorimetric	Green
	AS-71140	SensoLyte® Red Protease Assay Kit	Fluorimetric	Orange
Renin	AS-72040	SensoLyte® 520 Renin Activity Assay Kit	Fluorimetric	Green
	AS-72140	SensoLyte® 520 Rat Renin Activity Assay Kit	Fluorimetric	Green
	AS-72161	SensoLyte® 520 Mouse Renin Activity Assay Kit	Fluorimetric	Green
Sirtuin (SIRT)	AS-72155	SensoLyte® 520 FRET Sirt1 Activity Assay Kit	Fluorimetric	Green
	AS-72156	SensoLyte® Green Sirt1 Activity Assay Kit	Fluorimetric	Green
	AS-72189	SensoLyte® 520 FRET SIRT2 Activity Assay Kit	Fluorimetric	Green
	AS-72188	SensoLyte® Green SIRT2 Activity Assay Kit	Fluorimetric	Green
Sortase A	AS-72228	SensoLyte® 520 Sortase A Activity Assay Kit	Fluorimetric	Green
Thrombin	AS-72129	SensoLyte® 520 Thrombin Activity Assay Kit	Fluorimetric	Green
	AS-72130	SensoLyte® AFC Thrombin Activity Assay Kit	Fluorimetric	Blue
Tobacco Etch Virus (TEV) Protease	AS-72227	SensoLyte® 520 TEV Protease Activity Assay Kit	Fluorimetric	Green
Transglutaminase	AS-72244	SensoLyte® Transglutaminase Activity Assay Kit	Colorimetric	n/a
West Nile Virus (WNV) Protease	AS-72079	SensoLyte® 440 West Nile Virus Protease Activity Assay Kit	Fluorimetric	Blue
	AS-72080	SensoLyte® 570 West Nile Virus Protease Activity Assay Kit	Fluorimetric	Orange

## ELISA Kits

Target	Catalog #	Product name	Detection
Beta-Amyloid (1-40)	AS-55551	SensoLyte® Anti-Human $\beta$ -Amyloid (1-40) Quantitative ELISA Kit	Colorimetric
	AS-55553	SensoLyte® Anti-Mouse/Rat $\beta$ -Amyloid (1-40) Quantitative ELISA Kit	Colorimetric
Beta-Amyloid (1-42)	AS-55552	SensoLyte® Anti-Human $\beta$ -Amyloid (1-42) Quantitative ELISA Kit	Colorimetric
	AS-55554	SensoLyte® Anti-Mouse/Rat $\beta$ -Amyloid (1-42) Quantitative ELISA Kit	Colorimetric
Myelin Oligodendrocyte Glycoprotein (MOG)	AS-54465	SensoLyte® Anti-MOG(35-55) IgG Quantitative ELISA Kit (mouse/rat)	Colorimetric
	AS-55153-H	SensoLyte® Anti-Human MOG (1-125) Human IgG Specific ELISA Kit	Colorimetric
	AS-55153-M	SensoLyte® Anti-Human MOG (1-125) Mouse IgG Specific ELISA Kit	Colorimetric
	AS-55156	SensoLyte® Anti-Mouse MOG (1-125) IgG Quantitative ELISA Kit	Colorimetric
	AS-55157	SensoLyte® Anti-Rat MOG (1-125) IgG Quantitative ELISA Kit	Colorimetric
Myelin Proteolipid Protein (PLP)	AS-55523	SensoLyte® Anti-PLP (178-191) IgG Quantitative ELISA Kit (mouse)	Colorimetric
	AS-55524	SensoLyte® Anti-PLP (139-151) IgG Quantitative ELISA Kit (mouse)	Colorimetric
Synuclein	AS-55550-H	SensoLyte® Anti- $\alpha$ -Synuclein (Human) ELISA Kit	Colorimetric
	AS-55550-R	SensoLyte® Anti- $\alpha$ -Synuclein (Rat) ELISA Kit	Colorimetric

## Kits for ELISA and WB detection

Target	Catalog #	Product name	Detection	Color
Peroxidase	AS-71111	SensoLyte® ADHP Peroxidase Assay Kit	Fluorimetric	Orange
	AS-72127	SensoLyte® Luminescent Peroxidase Assay Kit	Luminometric	n/a
Alkaline phosphatase	AS-72146	SensoLyte® pNPP Alkaline Phosphatase Assay Kit	Colorimetric	n/a
	AS-71109	SensoLyte® FDP Alkaline Phosphatase Assay Kit	Fluorimetric	Green
	AS-72122	SensoLyte® Luminescent Alkaline Phosphatase Assay Kit	Luminometric	n/a

## Analyte Quantitation Kits

Target	Catalog #	Product name	Detection	Color
Biotin	AS-72096-120	SensoLyte® HABA Biotin Quantitation Kit	Colorimetric	n/a
	AS-72096-500	SensoLyte® HABA Biotin Quantitation Kit - Large size	Colorimetric	n/a
	AS-72163	SensoLyte® Biotin Quantitation Kit	Fluorimetric	Green
Glutathione	AS-72153	SensoLyte® Total GSH Assay Kit	Colorimetric	n/a
	AS-72154	SensoLyte® 520 Total GSH Assay kit	Fluorimetric	Green
	AS-72158	SensoLyte® Glutathione Cellular Assay Kit	Fluorimetric	Blue
Hydrogen Peroxide	AS-71112	SensoLyte® ADHP Hydrogen Peroxide Assay Kit	Fluorimetric	Orange
NADP/NADPH	AS-72205	SensoLyte® NADP/NADPH Assay Kit	Colorimetric	n/a
Phosphate	AS-71103	SensoLyte® MG Phosphate Assay Kit	Colorimetric	n/a
Thiol	AS-72136	SensoLyte® Thiol Quantitation Assay Kit	Colorimetric	n/a
	AS-72137	SensoLyte® ABD-F Thiol Quantitation Assay Kit	Fluorimetric	Blue
	AS-72138	SensoLyte® 520 Thiol Quantitation Assay Kit	Fluorimetric	Green

## Protein Quantitation Kit

Target	Catalog #	Product name	Detection	Color
Protein content	AS-71015	SensoLyte® OPA Protein Quantitation Kit	Fluorimetric	Blue

## Reporter Gene Assays

Target	Catalog #	Product name	Detection	Color
Beta-Galactosidase	AS-72132	SensoLyte® MUG b-Galactosidase Assay Kit	Fluorimetric	Blue
	AS-72133	SensoLyte® FDG b-Galactosidase Assay Kit	Fluorimetric	Green
	AS-72134	SensoLyte® ONPG b-Galactosidase Assay Kit	Colorimetric	n/a
Alkaline phosphatase	AS-72144	SensoLyte® pNPP Secreted Alkaline Phosphatase Reporter Gene Assay Kit	Colorimetric	n/a
	AS-72145	SensoLyte® Luminescent Secreted Alkaline Phosphatase Reporter Gene Assay Kit	Luminometric	n/a

## Amyloid Aggregation-related Products

Target	Catalog #	Product name	Detection	Color
Beta-Amyloid	AS-72213	SensoLyte® Thioflavin T β-Amyloid (1-40) Aggregation Kit	Fluorimetric	Blue
	AS-72214	SensoLyte® Thioflavin T β-Amyloid (1-42) Aggregation Kit	Fluorimetric	Blue
	AS-72215	AggreSure β-Amyloid (1-40) peptide	n/a	n/a
	AS-72216	AggreSure β-Amyloid (1-42) peptide	n/a	n/a
	AS-72071	SensoLyte Fluor Aβ(1-42) sampler kit	n/a	n/a