

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

## Enzyme Activity Assay kits

| Target                                         | Catalog # | Product name                                                                     | Detection    | Color  |
|------------------------------------------------|-----------|----------------------------------------------------------------------------------|--------------|--------|
| 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  |
| Meprin                                         | AS-72253  | Sensolyte® 520 Meprin α Activity Assay Kit                                       | Fluorimetric | Green  |
|                                                | AS-72254  | Sensolyte® 520 Meprin β Activity Assay Kit                                       | Fluorimetric | Green  |
| Neprilysin                                     | AS-72223  | Sensolyte® 520 Neprilysin 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  |

| Target                            | Catalog # | Product name                                               | Detection    | Color  |
|-----------------------------------|-----------|------------------------------------------------------------|--------------|--------|
| Peptidyl-prolyl isomerase (Pin1)  | AS-72240  | Sensolyte® Green PIN1 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 Fluo AB(1-42) sampler kit                      | n/a          | n/a   |