

Assay Kits



AnaSpec is proud to be the first in the industry to provide long wavelength FRET assays, HDAC assays; as well as anti-MOG (35-55) IgG Quantitative ELISA assays. Visit online for more detailed information of all our assay kits.

Anti-MOG (35-55) IgG Quantitative ELISA Assay Kits (mouse/rat)

| Detects | Total | Different IgG Subtypes | IgG ₁ | IgG _{2a} | IgG _{2b} | IgG _{2c} | IgG ₃ |
|---------|-------|------------------------|------------------|-------------------|-------------------|-------------------|------------------|
| Cat# | 54465 | 54471 | 54466 | 54467 | 54468 | 54469 | 54470 |

Cell Proliferation and Cytotoxicity Assay Kits

| Products | DHL™ Cell Cytotoxicity Assay Kit Ex/Em=560 nm/590 nm | DHL™ Cell Viability & Proliferation Assay Kit Ex/Em=560 nm/590 nm |
|--|---|--|
| Cell Proliferation & Cytotoxicity Assay Kits | 71302 (500 assays), 71303 (5000 assays) | 71300 (2000 assays), 71301 (10000 assays) |

HDAC Assay Kits

| Products | SensoLyte™ 440 HDAC Activity Assay Kit Ex/Em=354 nm/442 nm | SensoLyte™ 520 HDAC Activity Assay Kit Ex/Em=490 nm/520 nm |
|--------------------------|---|---|
| HDAC Activity Assay Kits | 72083 | 72084 |

Peroxidase Assay Kits

| Products | ADHP Hydrogen Peroxide Assay Kit Ex/Em=573 nm/590 nm | ADHP Peroxidase Assay Kit Ex/Em=573 nm/590 nm | ADHP Peroxidase ELISA Assay Kit Ex/Em=573 nm/590 nm |
|----------------------------|---|--|--|
| SensoLyte™ Peroxidase Kits | 71112 | 71111 | 71110-M, 71110-R |

Phosphatase Assay Kits

| Products | <i>Fluorimetric</i> | | <i>Colorimetric</i> |
|-------------------------------|---------------------|-------|---------------------|
| | FDP | MFP | pNPP |
| Kits | | | |
| Protein Phosphatase | 71100 | 71104 | 71105 |
| Alkaline Phosphatase | 71109 | | 71230 |
| Alkaline Phosphatase ELISA | 71101-M 71101-R | | 71232-M 71232-R |
| Secreted Alkaline Phosphatase | 71108 | | 71233 |
| Acid Phosphatase | | 71231 | |
| MG Phosphate | | | 71103 |

Protein Kinase Assay Kits

| | |
|---|--------------|
| Protein Kinase Substrate Sampler Kit * <i>Biotinylated</i> * | 71200 |
| Serine/Threonine Protein Kinase Substrate Sampler Kits * <i>FAM-labeled</i> * | 71201, 71202 |
| Serine/Threonine Protein Kinase Substrate Sampler Kits * <i>TAMRA-labeled</i> * | 72029, 72030 |
| Tyrosine Protein Kinase Substrate Profiling Kit * <i>Fluorimetric</i> * | 71203 |

Quantitation Assay Kits

| Products | AnaLyte™ OPA Protein Quantitation Kit Ex/Em=338 nm/455 nm | HABA Biotin Quantitation Kit * <i>Colorimetric</i> * Abs=500 nm |
|------------|--|--|
| Assay Kits | 71015 | 72096-120, 72096-500 |

Custom Service Available
 Assay Development
 Inhibitor Screening - HDAC, HIV, HCV, MMP & others

Protease Assay Kits

| | | | |
|------------|-------------|---|--|
| Cathepsins | Products | SensoLyte™ Cathepsin Assay Kits | SensoLyte™ 520 Cathepsin Assay Kits Ex/Em=490 nm/520 nm |
| | Cathepsin D | SensoLyte™ 390 Assay Kit (Ex/Em=330 nm/390 nm) 72098 | 72097 |
| | Cathepsin S | SensoLyte™ 440 Assay Kit (Ex/Em=354 nm/442 nm) 72100 | 72099 |

| | | | | | |
|---------|-------------------|--|--|---|---|
| Caspase | Products | SensoLyte™ AMC Series Ex/Em=340 nm/440 nm | SensoLyte™ AFC Series Ex/Em=370 nm/500 nm | SensoLyte™ Rh 110 Series Ex/Em=501 nm/527 nm | SensoLyte™ AnaRed Series Ex/Em=595 nm/635 nm |
| | Caspase 3/7 | 71118 | 71114 | 71141 | 71122 |
| | Caspase 8 | | 72088-100 (100 assays) 72088-200 (200 assays) | | |
| | Profiling | 71120 | 71116 | | |
| | Substrate Sampler | 71121 | 71117 | | |

| | | | | |
|---------|----------|---------------------------------------|---------------------------------------|---------------------------------------|
| HCV/HIV | Products | SensoLyte™ 490 Ex/Em=340 nm/440 nm | SensoLyte™ 520 Ex/Em=490 nm/520 nm | SensoLyte™ 620 Ex/Em=590 nm/620 nm |
| | HCV | 71126, 72087 | 71145 | 71146 |
| | HIV | 71127 | 71147 | 71148 |

| | | | | | |
|----------------|----------|-------------------------------------|--|--|---|
| MMP Assay Kits | Products | SensoLyte™ ELISA Kits Abs=450 nm | SensoLyte™ 490 Kits Ex/Em=340 nm/490 nm | SensoLyte™ 520 Kits Ex/Em=490 nm/520 nm | SensoLyte Plus™ 520 Kits Ex/Em=490 nm/520 nm |
| | MMP-1 | 72102 | 71128 | 71150 | 72012 |
| | MMP-2 | | 71129 | 71151 | |
| | MMP-3 | 72103 | 71130 | 71152 | |
| | MMP-7 | | 71132 | 71153 | |
| | MMP-8 | 72104 | 71133 | 71154 | |
| | MMP-9 | 72105 | 71134 | 71155 | 72017 |
| | MMP-10 | 72106 | | 72024 | |
| | MMP-12 | | 71137 | 71157 | |
| | MMP-13 | 72107 | 71135 | 71156 | 72019 |
| | MMP-14 | | | 72025 | |

SensoLyte™ 520 Generic MMP Activity Kit *Fluorimetric* - cat# 71158
 SensoLyte™ 570 Generic MMP Activity Kit *Fluorimetric* - cat# 72101
 SensoLyte™ Generic MMP Activity Kit *Colorimetric* - cat# 72095

SensoLyte™ 520 MMP Profiling Kit *Fluorimetric* - cat# 71136
 SensoLyte™ 520 MMP Substrate Sampler Kit *Fluorimetric* -
 cat# 71170

| | | |
|------------|----------------------------|---|
| Secretases | Products | SensoLyte™ 520 Assay Kit Ex/Em=490 nm/520 nm |
| | α-Secretase/TACE Assay Kit | 72085 |
| | β-Secretase Assay Kit | 71144 |

| | | | |
|-----|----------------------------------|---|---|
| WNV | Products | SensoLyte™ 440 Assay Kit Ex/Em=354 nm/442 nm | SensoLyte™ 570 Assay Kit Ex/Em=540 nm/575 nm |
| | West Nile Virus Assay Kits | 72079 | 72080 |

| | | | |
|-------------------|--|---|---|
| Renin-Angiotensin | Products | SensoLyte™ 390 Assay Kit Ex/Em=330 nm/390 nm | SensoLyte™ 520 Assay Kit Ex/Em=490 nm/520 nm |
| | Renin Assay Kit | 72039 | 72040 |
| | ACE2 (Angiotensin Converting Enzyme 2) Assay Kit | 72086 | |

| | | | |
|------------------|-----------------------------|--|--|
| Generic Protease | Products | SensoLyte™ Green Protease Assay Kit Ex/Em=488 nm/520 nm | SensoLyte™ Red Protease Assay Kit Ex/Em=546 nm/575 nm |
| | Generic Protease Assay Kits | 71124 | 71140 |