



integrated proteomics solutions.

PHOSPHATASE ASSAY KITS

AnaSpec, Inc. • 2149 O'Toole Ave. • San Jose, CA 95131 • 800-452-5530 • Fax 408-452-5059 • service@anaspec.com • www.anaspec.com

Still using *p*NPP to detect phosphatase activity? It is time to switch over to SensoLyte™ assays!

AnaSpec introduces a series of FDP and MFP-based fluorimetric assays to facilitate your research needs in phosphatase measurement. These fluorimetric assays are 100 times more sensitive than their colorimetric counterparts, and have a wider dynamic range. The fluorescence signal can easily be detected using the popular green fluorescence filter (Excitation/Emission=490±20 nm/520±20 nm), and are in a convenient mix and read format, ideal for high throughput screening of phosphatase activities or inhibitors or high volume ELISAs.

SensoLyte™ FDP Protein Phosphatase Assay Kit (cat# 71100):

Highly sensitive. Ideal for assaying protein phosphatases which require pH ≥ 6.5 for optimal activities.

SensoLyte™ MFP Protein Phosphatase Assay Kit (cat# 71104):

Highly sensitive. Ideal for assaying protein phosphatases which require pH < 6.5 for optimal activities.

SensoLyte™ FDP Alkaline Phosphatase ELISA Assay Kit (cat# 71101):

Highly sensitive. Alkaline phosphatase conjugated secondary antibody is included in this kit.

SensoLyte™ FDP Secreted Alkaline Phosphatase Reporter Gene Assay Kit (cat# 71108):

Highly sensitive. It is ideal for measuring the activity of placental alkaline phosphatase, a popular reporter gene product in medium, serum, or cell lysate.

SensoLyte™ FDP Alkaline Phosphatase Assay Kit (cat# 71109):

Highly sensitive. Ideal for measuring alkaline phosphatase activity in biological samples and for alkaline phosphatase-based ELISA assays (alkaline phosphatase conjugated secondary antibody is not included in this kit).

SensoLyte™ FDP Alkaline Phosphatase ELISA Assay Kit (cat# 71101):

Highly sensitive. Similar to cat# 71109, except alkaline phosphatase conjugated secondary antibody is included. Ideal for measuring alkaline phosphatase activity in alkaline phosphatase-based ELISA assays.

SensoLyte™ MFP Acid Phosphatase Assay Kit (cat# 71231):

Highly sensitive. Ideal for measuring acid phosphatase activity in biological samples. MFP maintains its strong fluorescence in acidic buffer. Useful for continuous kinetic measurement.

SensoLyte™ MG Phosphate Assay Kit (cat# 71103):

Colorimetric assay. Can be combined with the phosphorylated peptide-based protein phosphatase assay to measure the activity of protein phosphatase. Widely used for measuring organic and inorganic phosphate existing in a variety of solutions.

We also provide cost-effective colorimetric assay kits for experiments requiring medium sensitivity.

SensoLyte™ *p*NPP Protein Phosphatase Assay Kit (cat# 71105)

SensoLyte™ *p*NPP Alkaline Phosphatase ELISA Assay Kit (cat# 71232)

SensoLyte™ *p*NPP Alkaline Phosphatase Assay Kit (cat# 71230)

SensoLyte™ *p*NPP Secreted Alkaline Phosphatase Reporter Gene Assay Kit (cat# 71233)