

PEGylated protein ELISA kit

Validated, Complete Assay for the Sensitive Detection of a Diverse Set of PEGs in Biological Samples

The PEGylated Protein ELISA kit is a competitive assay specific to the backbone of poly(ethylene glycol) (PEG) and has been validated for use with a wide variety of PEG molecules including linear and branched forms as well as free and conjugated forms. As expected, the assay is more sensitive to higher molecular weight PEG molecules due to the repeating characteristics of the backbone. The assay uses a monoclonal antibody to PEG pre-coated on a microtiter plate. The antibody binds in a competitive manner the PEG or PEGylated protein in the sample or the PEG covalently linked to biotin. After incubation at room temperature the excess reagents are washed away and a streptavidin HRP conjugate is added to bind to the biotinylated PEG bound to the antibody. After a second wash to remove excess HRP substrate is added. An HRP-catalyzed reaction generates a blue color in the solution. The reaction is stopped with stop solution and the resulting yellow color is read at 450nm. The amount of signal is inversely proportional to the concentration of PEG.

Poly(ethylene glycol) (PEG) is a widely used polymer in drug delivery systems and is often directly conjugated to a drug therapeutic. PEGylation of drug therapeutics is desirable as it has been found to increase the retention time, reduce the immunogenicity and increase stability towards metabolic enzymes. The PEGylated Protein ELISA kit is applicable for drug development and pharmaceutical manufacturing applications including drug formulations, pharmacokinetics analysis, drug comparison, lead candidate identification, lot release criteria and in-process QC studies. At this time, there are no long-term studies revealing the fate of this polymer in the body. The need for toxicology studies that would determine the accumulation of the metabolized PEGylated therapeutics in various tissues has been identified.

Citations: 15

[View Online »](#)

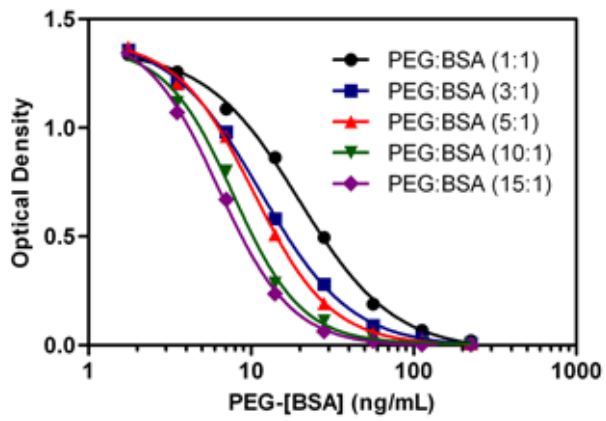
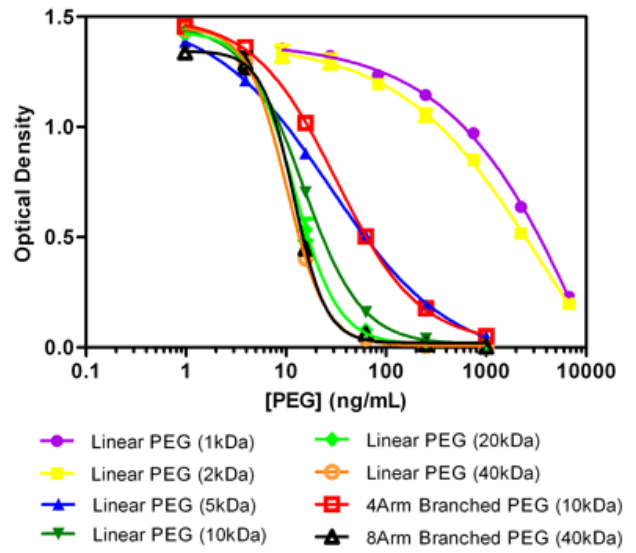
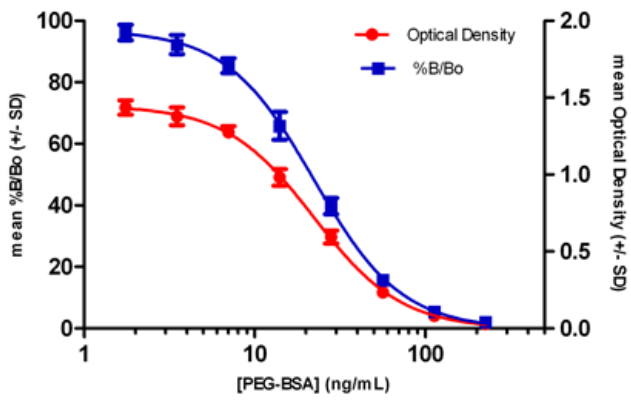
Ordering Information

[Order Online »](#)

ADI-900-213-0001

96 wells

- Validated for a wide range of MW linear and branched PEGs, both in free form and when conjugated to proteins
- Highly sensitive, measure as low as <1ng/mL of PEGylated proteins
- Quantify PEGylated target molecules in complex matrices to monitor drug levels or its accumulation in tissue
- High-throughput format, analyze up to 35 samples in duplicate results in just 2 hours



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Polyethylene glycol

Application Colorimetric detection, ELISA

Application Notes For the measurement of PEGylation and quantitative determination of PEGylated molecules in plasma, serum, tissue and other biological samples.

Assay Time 2 hours

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents Microtiter Plate, PEG-BSA Assay Control, Biotinylated PEG, Assay Buffer 39, Conjugate, Wash Buffer 3, TMB Substrate, Stop Solution 2, Plate Sealer

Species Reactivity Species independent

Wavelength 450 nm

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com