

Product datasheet

Polyethylene Glycol ELISA Kit ab215546

1 Image

Overview

Product name Polyethylene Glycol ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall	4			3%

Inter-assay

Sample	n	Mean	SD	CV%
Overall	4			6%

Sample type Serum, Plasma, Cell culture media

Assay type Competitive

Range 2 ng/ml - 200 ng/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	91	82% - 101%
Cell culture media	104	99% - 108%
Hep Plasma	70	64% - 77%
EDTA Plasma	84	77% - 90%
Cit plasma	79	75% - 85%

Assay duration Multiple steps standard assay

Product overview

Polyethylene Glycol (PEG) RabMAb[®] *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for accurate competitive quantitative measurement of PEGylated molecules in plasma, serum and cell culture media.

The Polyethylene Glycol (PEG) RabMAb[®] ELISA Kit operates on the basis of competition between enzyme HRP conjugated PEG and PEG labeled molecules for a limited number of binding sites on the surface of 96-wells coated with anti-PEG RabMAb[®] antibody. The extent of color development resulting from interaction between HRP and the substrate TMB is inversely proportional to the amount of PEGylated molecules in the sample. For example, the absence of PEGylated molecules in the sample will result in a bright blue color, whereas the presence of PEGylated molecules will result in decreased or no color development.

This kit is developed to determine levels of PEGylated molecules in samples such as serum, plasma or cell culture medium via ELISA.

Users must PEGylate their interested molecules. For pharmaco kinetic experiments, the PEGylated molecules will be used to construct standard curves.

The use of this kit requires the end user to have at least 250 ng of PEGylated compound of interest to use as reference standard. The included PEG-BSA is a reference sample only.

This kit does not contain Tween 20.

Consult our PEG FAQs page for more information:<https://www.abcam.com/primary-antibodies/polyethylene-glycol-peg-faqs>

Polyethylene Glycol ELISA Kit (ab138914) has been re-developed. Specifically, the way we manufacture some of the kit's components has changed (wash buffer, TMB and stop solution). This new kit has the same name but a different product number (ab215546), and the performance of the new kit is the same.

Notes Polyethylene glycol (PEG) is an O-CH₂-CH₂ polymer, which is water-soluble, nontoxic, non-antigenic, and biocompatible. Covalent conjugation of PEG to therapeutic proteins increases the *in vivo* stability by protecting the protein from degradation, masking its immunogenic sites and reducing clearance. Typically, PEGylation uses nonspecific reactions with nucleophilic residues and produces mixtures of PEGylated positional isomers. Qualitative and quantitative analysis of PEGylated molecules is important for both drug development and clinical application.

Platform Pre-coated microplate (12 x 8 well strips)

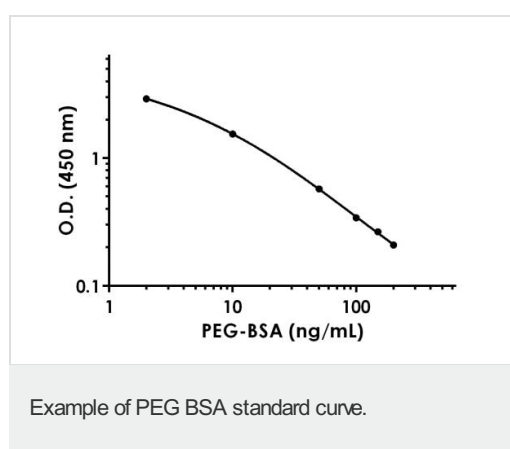
Properties

Storage instructions Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
10X Wash Buffer PT (ab206977)	1 x 20ml
Antigen/Antibody Diluent Buffer 1X	1 x 20ml
Anti-PEG coated microplate (12 x 8 well strips)	1 unit

Components	1 x 96 tests
PEG-conjugated HRP 60X	1 x 150µl
Plate Seals	1 unit
Reference Standard (PEG-BSA)	1 x 200ng
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

Images



Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors