

Mouse IGFBP-1 ELISA Kit ab213865

1 References 1 Image

Overview

Product name	Mouse IGFBP-1 ELISA Kit			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	1	16	88pg/ml	6.6
	2	16	281pg/ml	14.77
	3	16	865pg/ml	38.06
	Inter-assay			
	Sample	n	Mean	SD
	1	24	90pg/ml	7.92
	2	24	267pg/ml	19.49
	3	24	783pg/ml	36.8
Sample type	Cell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma, Tissue Homogenate			
Assay type	Sandwich (quantitative)			
Sensitivity	< 3 pg/ml			
Range	31.2 pg/ml - 2000 pg/ml			
Assay time	3h 30m			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Mouse			
Product overview	The Mouse IGFBP-1 Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213865) is designed for the quantitative detection of Mouse IGFBP-1 in cell culture supernatants, cell lysates, tissue homogenates, serum and plasma (heparin, EDTA).			

The ELISA kit is based on standard sandwich enzyme-linked immunosorbent assay technology. A

polyclonal antibody from goat specific for IGFBP-1 has been pre-coated onto 96-well plates. Standards (Expression system for standard: NSO; Immunogen sequence: A26-N272) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for IGFBP-1 is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex was added and unbound conjugates were washed away with PBS or TBS buffer. HRP substrate TMB was used to visualize HRP enzymatic reaction. TMB was catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the Mouse IGFBP-1 amount of sample captured in plate. The standard product used in this kit is recombinant mouse IGFBP-1, consisting of 247 amino acids with the molecular mass of 27 KDa.

Notes IGFBP-1 is also known as amniotic fluid binding protein (AFBP), placental protein-12, alpha-pregnancy associated endometrial globulin, growth hormone independent binding protein, binding protein-28, binding protein-26, and binding protein-25. The low-molecular weight insulin-like growth-factor binding protein (IGF-BP25) is synthesized by human liver, secretory endometrium and decidua, and is also present in human serum. It binds insulin-like growth factors IGF-I and IGF-II with high affinity, and is proposed to act as a paracrine regulator of cell growth. These IGF-binding proteins are expressed at different concentrations in different tissues and are thought to regulate the activity of IGF I and II. The gene is organized in four exons and spans 5.9 kb. The IBP-1 gene is a single copy gene, located on chromosome 7.

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

Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	Identifier	1 x 96 tests
TMB Stop Solution		—
Avidin-Biotin-Peroxidase Complex (ABC)		—
Sample Diluent Buffer		—
Antibody Diluent Buffer		—
ABC Diluent Buffer		—
TMB Color Developing Agent		—
Adhesive Plate Seal		—
Anti-Mouse IGFBP-1 coated Microplate (12 x 8 wells)		—
Lyophilized recombinant Mouse IGFBP-1 standard		—
Biotinylated anti- Mouse IGFBP-1 antibody		—
Wash Buffer (25X)		—

Function IGF-binding proteins prolong the half-life of the IGFs and have been shown to either inhibit or stimulate the growth promoting effects of the IGFs on cell culture. They alter the interaction of IGFs

with their cell surface receptors. Promotes cell migration.

Sequence similarities

Contains 1 IGFBP N-terminal domain.

Contains 1 thyroglobulin type-1 domain.

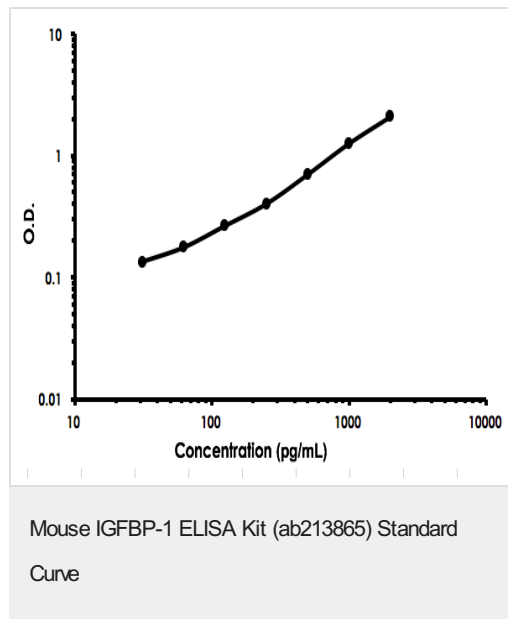
Post-translational modifications

Phosphorylated; probably by casein kinase II. Alters the affinity of the protein for IGFs.

Cellular localization

Secreted.

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