

Product datasheet

Mouse IGFBP6 Matched Antibody Pair Kit ab215412

1 Image

Overview

Product name	Mouse IGFBP6 Matched Antibody Pair Kit
Detection method	Colorimetric
Assay type	ELISA set
Sensitivity	4.4 pg/ml
Range	31.25 pg/ml - 2000 pg/ml
Species reactivity	Reacts with: Mouse
Product overview	<p>Mouse IGFBP6 Matched Antibody Pair Kits include a capture and a biotinylated detector antibody pair, along with a calibrated protein standard, suitable for sandwich ELISA. The Matched Antibody Pair Kit can be used to quantify native and recombinant mouse IGFBP6.</p>

Optimization of the kit reagents to sample type, immunoassay format or instrumentation may be required. Guidelines for use of this kit in a standard 96-well microplate sandwich ELISA using HRP/TMB system of colorimetric detection is described in this assay procedure for the purposes of quantification.

Protocol information and tips on the use of the Matched Antibody Pair kits for sandwich ELISA can be found on our [website](#). An accessory pack can be purchased which includes buffer reagents required to perform 10 x 96-well plate sandwich ELISAs ([ab210905](#)).

For additional information on the performance of the antibody pair used in this kit, please see our equivalent SimpleStep ELISA kit [ab206305](#). Please note that while the antibody pair is the same provided in the corresponding SimpleStep ELISA Kit, due to differences in their formulation, this antibody pair cannot be used with the consumables provided with our SimpleStep ELISA Kits.

We've listened to you: due to popular demand, we will now provide our Matched Antibody Pair kits in 5x96 tests and 10x96 tests. The 2x96 tests size will be discontinued on 30th June 2020 unless inventory is depleted beforehand.

To receive an electronic copy of the Certificate of Analysis, please send an [email](#) with "CoA for matched antibody pair kit" in the subject line and the desired product number and lot number in the body of the email.

Tested applications	Suitable for: IA, ELISA
Platform	Reagents

Properties

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

Components	10 x 96 tests	5 x 96 tests
Mouse IGFBP6 Capture Antibody	1 x 100µg	1 x 50µg
Mouse IGFBP6 Detector Antibody	1 x 25µg	1 x 12.5µg
Mouse IGFBP6 Lyophilized Protein	1 vial	1 vial

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.

Sequence similarities Contains 1 IGFBP N-terminal domain.
Contains 1 thyroglobulin type-1 domain.

Post-translational modifications O-linked glycans consist of hexose (probably Gal), N-acetylhexosamine (probably GalNAc) and sialic acid residues. O-glycosylated with core 1 or possibly core 8 glycans. O-glycosylated on one site only in the region AA 143-168 in cerebrospinal fluid.

Cellular localization Secreted.

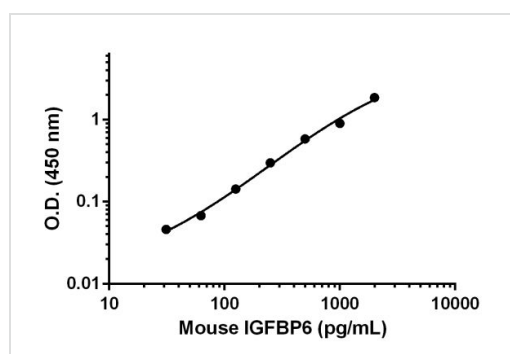
Applications

Our [Abpromise guarantee](#) covers the use of **ab215412** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
IA		Use at an assay dependent concentration.
ELISA		Use at an assay dependent concentration.

Images



Mouse IGFBP6 standard curve.

Standard calibration curve. Background subtracted values are graphed.

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