

Product datasheet

Mouse IGFBP2 Matched Antibody Pair Kit ab216787

1 Image

Overview

| | |
|---------------------------|--|
| Product name | Mouse IGFBP2 Matched Antibody Pair Kit |
| Detection method | Colorimetric |
| Assay type | ELISA set |
| Range | 93.75 pg/ml - 6000 pg/ml |
| Species reactivity | Reacts with: Mouse |
| Product overview | Matched Antibody Pair kit is a titrated unlabeled capture antibody, a titrated biotin-labeled detector and a calibrated protein standard. The Matched Antibody Pair Kit can be used to quantify native and recombinant mouse IGFBP2. |

Both capture and detector antibodies are rabbit monoclonal antibodies.

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.

Additional protocol information and tips on the use of the Matched Antibody Pair kits for sandwich ELISA can be found on our [website](#).

For additional information on the performance of the antibody pair used in this kit, please see our equivalent SimpleStep ELISA® ([ab207615](#)) which uses the same antibody pair.

Buffer information: The capture antibody is glycerol free. The detector antibody contains glycerol.

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: ELISA |
| Platform | Reagents |

Properties

Storage instructions

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

| Components | 10 x 96 tests | 2 x 96 tests |
|----------------------------------|---------------|--------------|
| Mouse IGFBP2 Capture Antibody | 1 x 100µg | 1 x 20µg |
| Mouse IGFBP2 Detector Antibody | 1 x 25µg | 1 x 5µg |
| Mouse IGFBP2 Lyophilized Protein | 1 vial | 1 vial |

Function

Inhibits IGF-mediated growth and developmental rates. 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.

Domain

The C-terminus is required for IGF-binding and growth inhibition.

Post-translational modifications

O-glycosylated.

Cellular localization

Secreted.

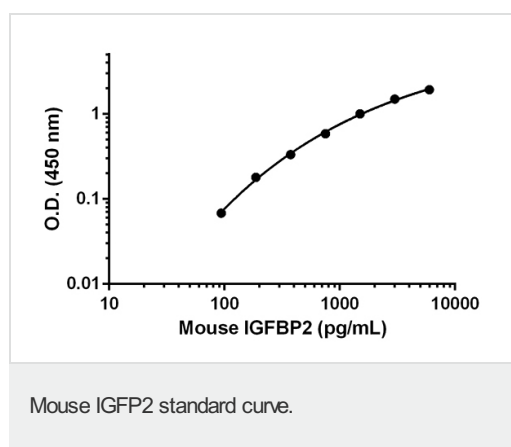
Applications

Our [Abpromise guarantee](#) covers the use of **ab216787** 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 |
|-------------|-----------|--|
| ELISA | | Use at an assay dependent concentration. |

Images



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