

Mouse IGFBP2 ELISA Kit ab100691

[1 References](#) [1 Image](#)

Overview

Product name	Mouse IGFBP2 ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 0.6 ng/ml
Range	0.823 ng/ml - 600 ng/ml
Recovery	91 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	112.5	103% - 122%
Serum	81.45	72% - 90%
Plasma	80.22	73% - 92%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview Abcam's IGFBP2 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Mouse IGFBP2 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Mouse IGFBP2 coated on a 96- well plate. Standards and samples are pipetted into the wells and IGFBP2 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Mouse IGFBP2 antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of IGFBP2 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Platform Microplate

Properties

Storage instructions

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

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
500X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent C	1 x 30ml
Biotinylated anti-Mouse IGFBP2	2 vials
IGFBP2 Microplate (12 x 8 wells)	1 unit
Recombinant Mouse IGFBP2 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

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.

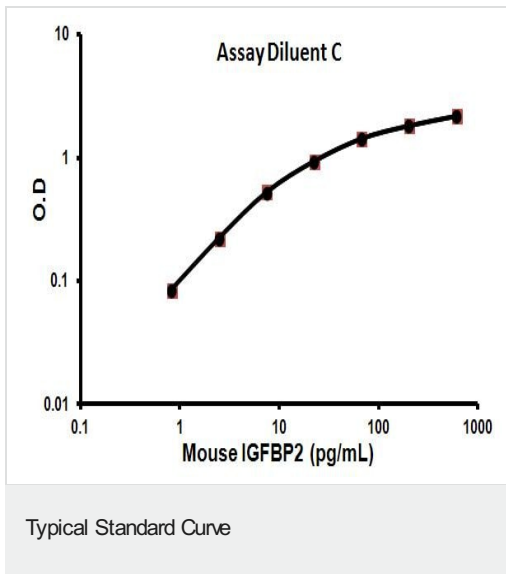
Post-translational modifications

O-glycosylated.

Cellular localization

Secreted.

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