

Mouse IGFBP5 ELISA Kit ab100693

1 References 1 Image

Overview

Product name	Mouse IGFBP5 ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 5 pg/ml
Range	4.12 pg/ml - 3000 pg/ml
Recovery	97 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	91.26	80% - 101%
Serum	115.3	86% - 130%
Plasma	95.11	83% - 102%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview Abcam's IGFBP5 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Mouse IGFBP5 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for mouse IGFBP5 coated on a 96- well plate. Standards and samples are pipetted into the wells and IGFBP5 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse IGFBP5 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 IGFBP5 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Get better reproducibility in only 90 minutes with Mouse IGFBP5 ELISA Kit ([ab208345](#)) from our

SimpleStep ELISA® range.

Platform

Microplate

Properties

Storage instructions

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

Components	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent C	1 x 30ml
Biotinylated anti-Mouse IGFBP5	2 vials
IGFBP5 Microplate (12 x 8 wells)	1 unit
Recombinant Mouse IGFBP5 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

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.

Tissue specificity

Osteosarcoma, and at lower levels in liver, kidney and brain.

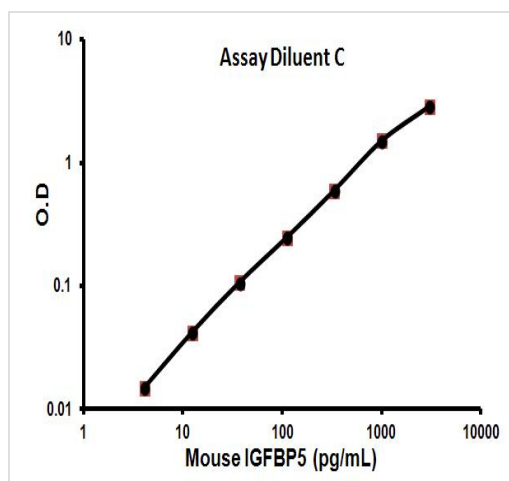
Sequence similarities

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

Cellular localization

Secreted.

Images



Representative Standard Curve using ab100693

Typical Standard Curve

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