

Mouse IGFBP2 ELISA Kit ab100691

[1 References](#) [1 图像](#)

概述

| | |
|------|---|
| 产品名称 | 小鼠IGFBP2 ELISA试剂盒 |
| 检测方法 | Colorimetric |
| 样品类型 | Cell culture supernatant, Serum, Plasma |
| 检测类型 | Sandwich (quantitative) |
| 灵敏度 | < 0.6 ng/ml |
| 范围 | 0.823 ng/ml - 600 ng/ml |
| 回收率 | 91 % |

特定样本回收率

| 样品类型 | 平均% | 范围 |
|--------------------------|-------|-------------|
| Cell culture supernatant | 112.5 | 103% - 122% |
| Serum | 81.45 | 72% - 90% |
| Plasma | 80.22 | 73% - 92% |

实验步骤 Multiple steps standard assay

种属反应性 与反应: Mouse

产品概述 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.

平台 Microplate

性能

存放说明

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

| 组件 | 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 |

功能

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.

序列相似性

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

结构域

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

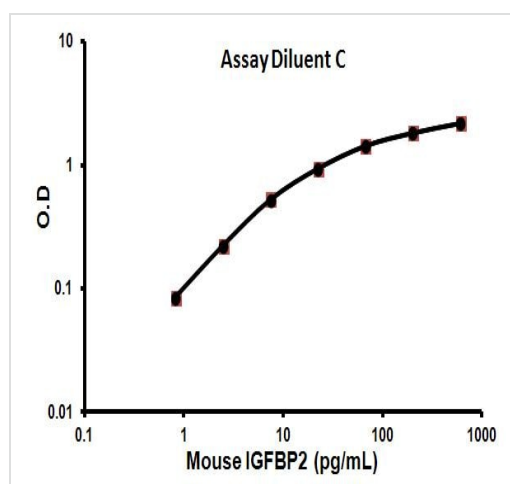
翻译后修饰

O-glycosylated.

细胞定位

Secreted.

图片



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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors