

Human IGFBP2 ELISA Kit ab100540

4 References 2 图像

概述

产品名称	人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	< 20 pg/ml
范围	8.23 pg/ml - 6000 pg/ml
回收率	97 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	100.25	86% - 105%
Serum	95.77	84% - 103%
Plasma	97.59	85% - 104%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

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

This assay employs an antibody specific for Human 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-Human 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.

说明 Optimization may be required with urine samples

平台 Microplate

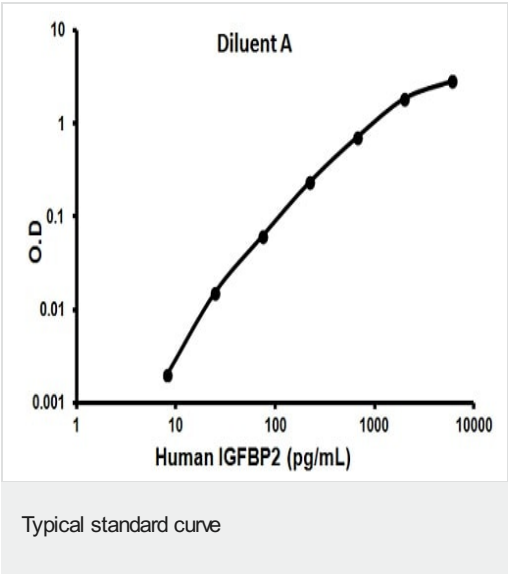
性能

存放说明Store at -20°C. Please refer to protocols.

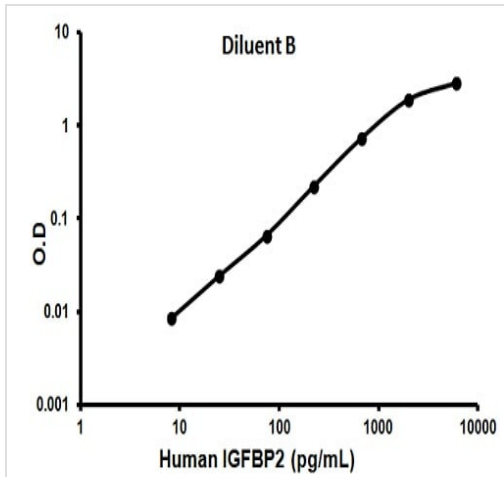
组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
300X HRP-Streptavidin Concentrate	1 x 200μl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human IGFBP2	2 vials
IGFBP2 Microplate (12 x 8 wells)	1 unit
Recombinant Human IGFBP2 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step 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



Representative standard curve using ab100540



Representative standard curve using ab100540

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