

Product datasheet

Human IGFBP3 ELISA Kit ab100541

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

Overview

Product name	Human IGFBP3 ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 80 pg/ml
Range	74.07 pg/ml - 18000 pg/ml
Recovery	97 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	97.46	86% - 107%
Serum	97.46	85% - 105%
Plasma	96.69	84% - 106%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's IGFBP3 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human IGFBP3 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Human IGFBP3 coated on a 96-well plate. Standards and samples are pipetted into the wells and IGFBP3 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human IGFBP3 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 IGFBP3 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Notes Optimization may be required with urine samples

Platform Microplate

Properties

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

Components	1 x 96 tests
20X Wash Buffer Concentrate	1 x 25ml
320X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human IGFBP3 (lyophilized)	2 vials
IGFBP3 Microplate (12 x 8 wells)	1 unit
Recombinant Human IGFBP3 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

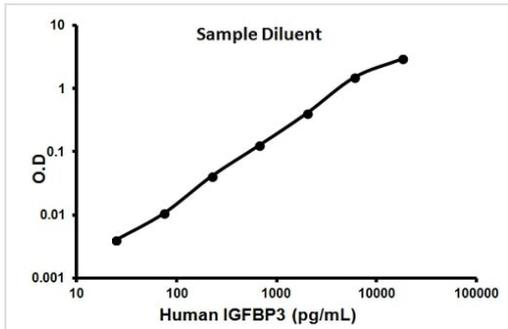
Relevance

Insulin like growth factor binding protein 3 (IGFBP3) is a member of the superfamily of insulin like growth factor (IGF) binding proteins which include six high affinity IGF binding proteins (IGFBP) and at least four low affinity binding proteins referred to as IGFBP related proteins (IGFBPrP). The IGFBP members are cysteine rich proteins with conserved cysteine residues clustered in the amino terminal and the carboxy terminal regions of the molecule. The cDNA sequence encoding the mature human IGFBP3 is fused to the signal peptide of CD33. Human IGFBP3 is the major IGF binding protein in plasma where it exists in a ternary complex with IGF1 or IGFII and an acid labile subunit. IGFBPs hold a central position in IGF ligand-receptor interactions through influences on both the bioavailability and distribution of IGFs in the extracellular environment. Insulin like growth factor binding protein 3 (IGFBP3) can modulate the mitogenic and metabolic effects of the insulin like growth factors (IGFs). Insulin like growth factor binding protein 3 (IGFBP3) is expressed in multiple tissues. The highest expression level is found in the non paranchymal cells of the liver. The expression levels are higher during extrauterine life and peak during puberty. Insulin like growth factors (IGFs) and IGF binding proteins (IGFBPs) play important roles in cell growth and differentiation. IGFBP3 is one of the factors in serum that is responsible for high serum induced apoptosis in PC3 cells, a prostate cancer cell line. IGFBP3 is important in controlling glucose homeostasis with increased urinary levels in type I diabetes with persistent microalbuminuria.

Cellular localization

Secreted

Images



Representative standard curve using ab100541

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