

Human IL-18BP ELISA Kit ab100559

3 References 2 Images

Overview

Product name	Human IL-18BP ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 20 pg/ml
Range	24.69 pg/ml - 18000 pg/ml
Recovery	91 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	106.7	96% - 121%
Serum	83.5	74% - 97%
Plasma	82.3	72% - 96%

Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Human
Product overview	Abcam's IL-18BP (Interleukin-18 Binding Protein) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human IL-18BP in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human IL-18BP coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-18BP present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human IL-18 BP 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 IL-18BP 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	1 x 25ml
5X Assay Diluent B	1 x 15ml
500X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated anti-Human IL-18BP	2 vials
IL-18BP Microplate (12 x 8 wells)	1 unit
Recombinant Human IL-18BP Standard (lyophilized)	2 units
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Function

Isoform A binds to IL-18 and inhibits its activity. Functions as an inhibitor of the early TH1 cytokine response.

Tissue specificity

Strongly expressed in heart, lung, placenta and spleen.

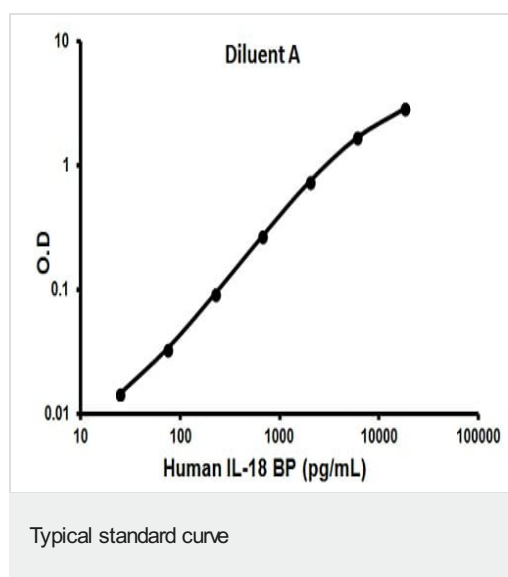
Sequence similarities

Contains 1 Ig-like C2-type (immunoglobulin-like) domain.

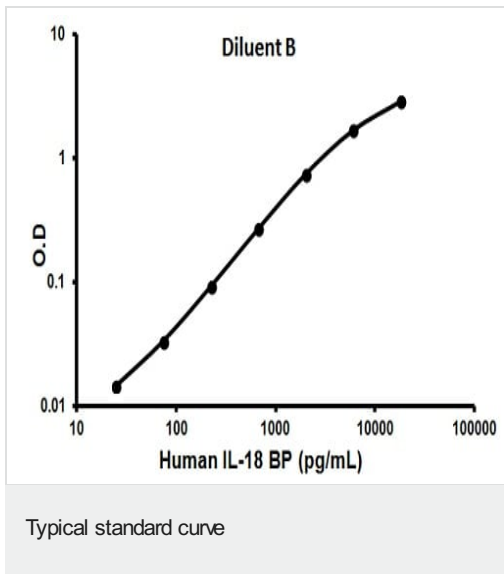
Cellular localization

Secreted.

Images



Representative standard curve using ab100559



Representative standard curve using ab100559

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