

Human IL-18BP ELISA Kit ab100559

[3 References](#) [2 图像](#)

概述

产品名称	人IL-18BP 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	24.69 pg/ml - 18000 pg/ml
回收率	91 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	106.7	96% - 121%
Serum	83.5	74% - 97%
Plasma	82.3	72% - 96%

实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	<p>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.</p> <p>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.</p>
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
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

功能

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

组织特异性

Strongly expressed in heart, lung, placenta and spleen.

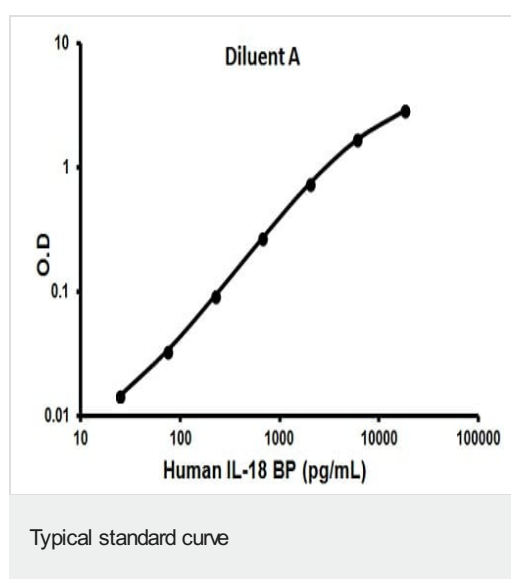
序列相似性

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

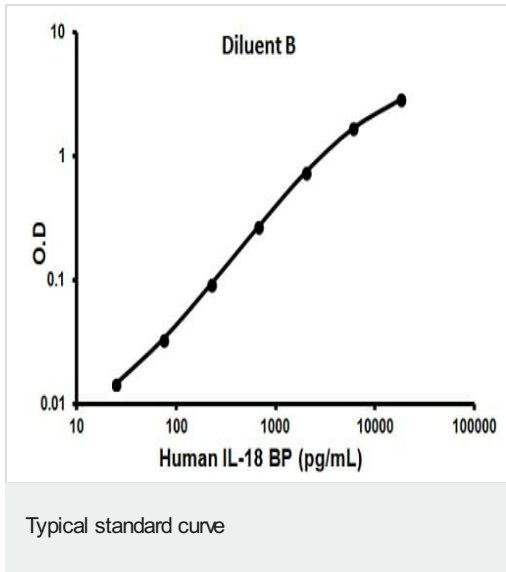
细胞定位

Secreted.

图片



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.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