

Human IL-18BP ELISA Kit ab100559

3 References 画像数 2

製品の概要

製品名	Human IL-18BP ELISA Kit
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 20 pg/ml
検出範囲	24.69 pg/ml - 18000 pg/ml
添加回収試験	91 %

特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	106.7	96% - 121%
Serum	83.5	74% - 97%
Plasma	82.3	72% - 96%

ステップ	Multiple steps standard assay
種交差性	交差種: Human
製品の概要	<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>
特記事項	Optimization may be required with urine samples.
試験プラットフォーム	Microplate

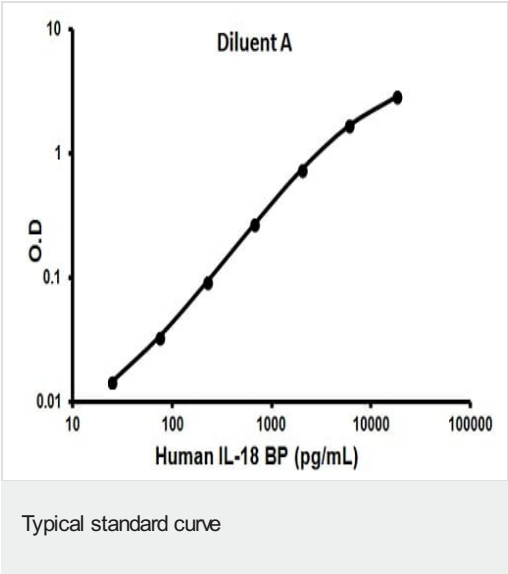
製品の特性

保存方法Store at -20°C. Please refer to protocols.

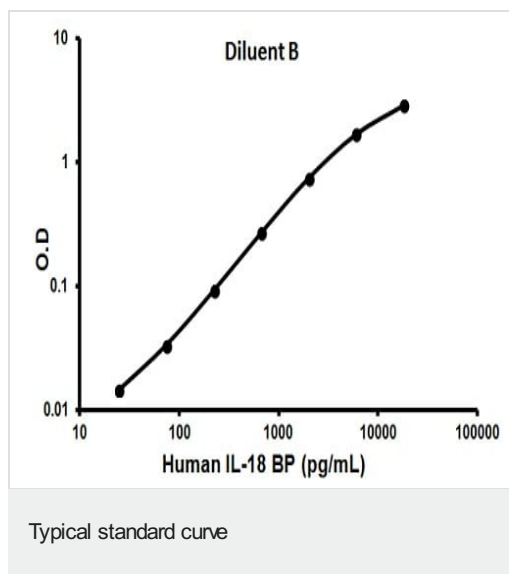
内容	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
500X HRP-Streptavidin Concentrate	1 x 200μl
Assay Diluent A	1 x 30ml
Biotinylated anti-Human IL-18BP	2 vials
IL-18BP Microplate (12 x 8 wells)	1 unit
Recombinant Human IL-18BP Standard (lyophilized)	2 units
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

機能	Isoform A binds to IL-18 and inhibits its activity. Functions as an inhibitor of the early TH1 cytokine response.
組織特異性	Strongly expressed in heart, lung, placenta and spleen.
配列類似性	Contains 1 Ig-like C2-type (immunoglobulin-like) domain.
細胞内局在	Secreted.

画像



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.co.jp/abpromise> or contact our technical team.

#### Terms and conditions

---

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