

Human IL-13 ELISA Kit ab100553

3 References 画像数 2

製品の概要

製品名	Human IL-13 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	< 0.15 pg/ml
検出範囲	0.16 pg/ml - 40 pg/ml
添加回収試験	100 %

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	100.65	88% - 107%
Serum	96.35	84% - 103%
Plasma	94.39	82% - 102%

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 Abcam's IL-13 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human IL-13 in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human IL-13 coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-13 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human IL-13 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-13 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Get results in 90 minutes with Human IL-13 ELISA Kit ([ab178014](#)) from our SimpleStep ELISA® range.

特記事項	Optimization may be required with urine samples
試験プラットフォーム	Microplate

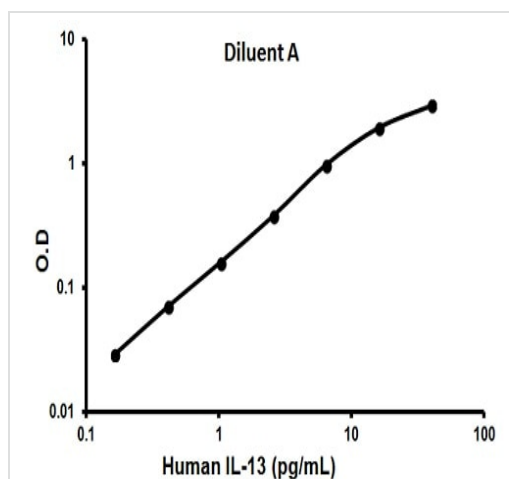
製品の特性

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

内 容	1 x 96 tests
20X Wash Buffer	1 x 25ml
120X HRP-Streptavidin Concentrate	1 x 200μl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human IL-13	2 vials
IL-13 Microplate (12 x 8 wells)	1 unit
Recombinant Human IL-13 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

機 能	Cytokine (PubMed:8096327, PubMed:8097324). Inhibits inflammatory cytokine production (PubMed:8096327). Synergizes with IL2 in regulating interferon-gamma synthesis (PubMed:8096327). May be critical in regulating inflammatory and immune responses (PubMed:8096327, PubMed:8097324). Positively regulates IL31RA expression in macrophages.
関連疾患	Allergic rhinitis
配列類似性	Belongs to the IL-4/IL-13 family.
細胞内局在	Secreted.

画 像



Representative standard curve using ab100553

Typical standard curve



Representative standard curve using ab100553

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.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