

Human IL-4 ELISA Kit ab46058

画像数 1

製品の概要

製品名 Human IL-4 ELISA Kit

検出方法 Colorimetric

再現性 Intra-Assay (同時再現性)

| サンプル | N | 平均値 | SD | CV% |
|--------|---|-----|----|------|
| sample | 6 | | | 3.5% |

Inter-Assay (日差再現性)

| サンプル | N | 平均値 | SD | CV% |
|--------|---|-----|----|------|
| sample | 6 | | | 7.3% |

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 < 0.7 pg/ml

検出範囲 1.1 pg/ml - 35 pg/ml

添加回収試験 107 %

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 |
|---------|------|------------|
| Serum | 100 | 75% - 121% |

全工程の試験時間 3h 45m

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 Abcam's IL-4 Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-4 in supernatants, buffered solutions, serum and plasma samples.

A monoclonal antibody specific IL-4 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IL-4 concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and

a biotinylated monoclonal antibody specific for IL-4 are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated and washed. A TMB substrate solution is added which acts on the bound enzyme to induce a colored reaction product. The intensity of this colored product is directly proportional to the concentration of IL4 present in the samples.

This kit will recognize both endogenous and recombinant Human IL-4.

試験プラットフォーム

Microplate

製品の特性

保存方法

Store at +4°C. Please refer to protocols.

| 内容 | ラベル | 1 x 96 tests | 2 x 96 tests |
|--------------------------------------|-------|--------------|--------------|
| 10X Standard Diluent Buffer | Black | 1 x 15ml | 1 x 25ml |
| 200X Wash Buffer | White | 1 x 10ml | 2 x 10ml |
| Biotinylated anti IL-4 | | 1 x 400µl | 2 x 400µl |
| Biotinylated Antibody Diluent | Red | 1 x 7ml | 1 x 13ml |
| Chromogen TMB Substrate Solution | | 1 x 11ml | 1 x 24ml |
| Control | | 2 vials | 4 vials |
| HRP Diluent | Red | 1 x 12ml | 1 x 23ml |
| IL-4 Microplate (12 x 8 well strips) | | 1 unit | 2 units |
| IL-4 Standard (Lyophilized) | | 2 vials | 4 vials |
| Standard Diluent (Serum) | | 1 x 7ml | 2 x 7ml |
| Stop Reagent | Black | 1 x 11ml | 2 x 11ml |
| Streptavidin-HRP | | 2 x 5µl | 4 x 5µl |

関連性

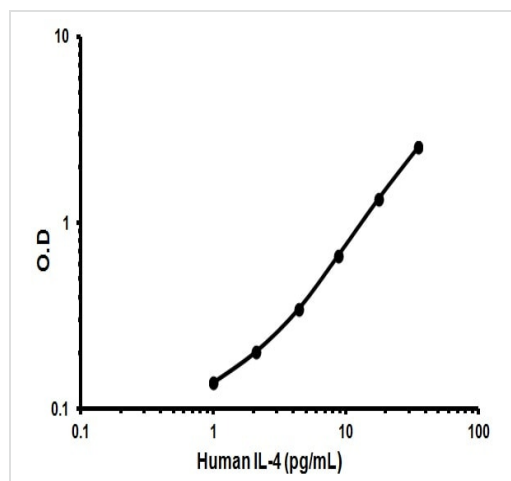
IL-4 is a pleiotropic cytokine produced by activated T cells. It is a ligand for interleukin 4 receptor. The interleukin 4 receptor also binds to IL13, which may contribute to many overlapping functions of this cytokine and IL13. STAT6, a signal transducer and activator of transcription, has been shown to play a central role in mediating the immune regulatory signal of IL-4. IL-4, IL3, IL5, IL13, and CSF2 form a cytokine gene cluster on chromosome 5q, with this gene particularly close to IL13. IL-4, IL13 and IL5 are found to be regulated coordinately by several long-range regulatory elements in an over 120 kilobase range on the chromosome. Two alternatively spliced transcript variants of IL-4 encoding distinct isoforms have been reported.

細胞内局在

Secreted

画像

Representative standard curve using ab46058



Typical data

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