

Human IL-4 ELISA Kit High Sensitivity ab46063

2 References [画像数 3](#)

製品の概要

製品名 Human IL-4 ELISA Kit High Sensitivity

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%
A	8	7.43pg/ml	0.9	12.3%
B	8	6.2pg/ml	0.7	10.8%
C	8	8.77pg/ml	0.7	7.6%
D	8	7.36pg/ml	0.4	6%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
A	9	7.1pg/ml	0.83	11.8%
B	9	6.47pg/ml	0.72	11.1%
C	9	8.76pg/ml	1.37	15.7%
D	9	7.44pg/ml	1.2	16.1%
E	7	2pg/ml	0.874	17%

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 = 0.31 pg/ml

検出範囲 0.31 pg/ml - 10 pg/ml

添加回収試験 101 %

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	95	97% - 127%
Serum	69	61% - 75%

全工程の試験時間

3h 00m

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 High Sensitivity *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-4 in Human serum, plasma, buffered solution and cell culture media.

A monoclonal antibody specific for 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 IL-4 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.

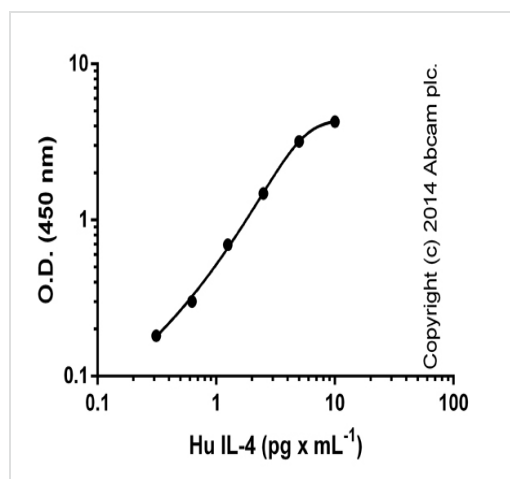
内容	ラベル	1 x 96 tests
10X Standard Diluent Buffer	Black	1 x 15ml
200X Wash Buffer	White	1 x 10ml
Amplification Diluent	Brown & blue spot	1 x 15ml
Amplifier	Yellow	1 x 200µl
Biotinylated anti IL-4		1 x 400µl
Biotinylated Antibody Diluent	Red	1 x 7.5ml
Chromogen TMB Substrate Solution		1 x 11ml
HRP Diluent	Red	1 x 25ml

内容	ラベル	1 x 96 tests
IL-4 Microplate (12 x 8 well strips)		1 unit
IL-4 Standard (Lyophilized)		2 vials
Stop Reagent	Black	1 x 11ml
Streptavidin-HRP		2 x 5 μ l

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.

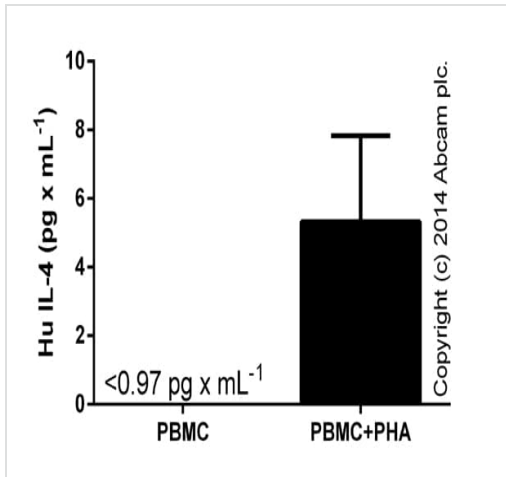
細胞内局在 Secreted

画像



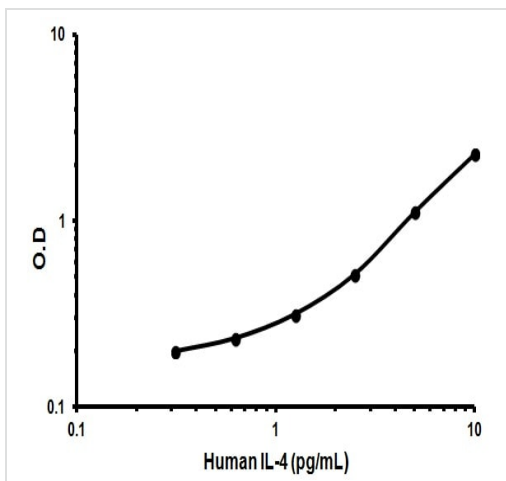
Standard curve of human IL-4, with background signal subtracted (duplicates; +/- SD).

ELISA: IL-4 High Sensitivity Human ELISA Kit
(ab46063)



IL-4 measured in cell culture supernatants from control (PBMC) or treated (PBMC+PHA; 2% PHA-M; LifeTechnologies) cells, grown for 48 hours (duplicates +/- SD). Samples tested neat - 1/10, control supernatant gave signals below level of detection.

ELISA: IL-4 High Sensitivity Human ELISA Kit (ab46063)



Representative Standard Curve using ab46063

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