

Product datasheet

Human IL-12 p70 ELISA Kit (Interleukin-12 p70) ab46143

画像数 1

製品の概要

製品名 Human IL-12 p70 ELISA Kit (Interleukin-12 p70)

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%
Overall				4.3%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Overall	6			8.9%

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 = 2.2 pg/ml

検出範囲 6.25 pg/ml - 200 pg/ml

添加回収試験 95 %

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	91	86% - 94%

全工程の試験時間 3h 40m

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要

Abcam's IL-12 p70 (Interleukin- 12 p70) Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-12 p70 in supernatants, buffered solutions or serum and plasma samples.

A monoclonal antibody specific IL-12 p70 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IL-12 p70 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-12 p70 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-12 p70 present in the samples.

This kit will recognize both endogenous and recombinant Human IL-12 p70.

アプリケーション

適用あり: Sandwich ELISA

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	2 x 96 tests	1 x 96 tests
10X Standard Diluent Buffer	Black	1 x 25ml	1 x 25ml
Wash Buffer	White	2 x 10ml	1 x 10ml
Biotinylated Antibody Diluent	Red	1 x 13ml	1 x 7.5ml
Biotinylated anti-IL-12 p70	Red	2 x 400µl	1 x 400µl
TMB Substrate		1 x 24ml	1 x 11ml
Control	Silver	4 vials	2 vials
HRP Diluent	Red	1 x 23ml	1 x 23ml
IL-12 p70 Microplate (12 x 8 well strips)		2 units	1 unit
IL12p70 Standard (Lyophilized)	Yellow	4 vials	2 vials
Standard Diluent (Human Serum)		1 x 7ml	2 x 7ml
Stop Reagent	Black	2 x 11ml	1 x 11ml
Streptavidin-HRP		4 x 5µl	2 x 5µl

機能

Cytokine that can act as a growth factor for activated T and NK cells, enhance the lytic activity of NK/lymphokine-activated Killer cells, and stimulate the production of IFN-gamma by resting PBMC.

配列類似性

Belongs to the IL-6 superfamily.

細胞内局在

Secreted.

アプリケーション

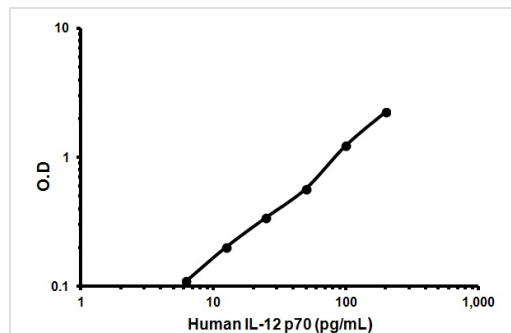
Our [Abpromise guarantee](#) covers the use of **ab46143** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Sandwich ELISA

Use at an assay dependent concentration.

画像



Typical Standard Curve

Representative Standard Curve using
ab46143

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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