

Human IL-12 ELISA Kit High Sensitivity ab46035

[5 References](#) [1 图像](#)

概述

产品名称 人IL-12 ELISA试剂盒High Sensitivity

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
Sample A	8	16.1pg/ml	1.25	7.8%
Sample B	8	6.1pg/ml	0.4	6.6%

批次间

样品	n	Mean	SD	CV%
Sample A	20	16pg/ml	1.2	7.9%
Sample B	20	6pg/ml	0.6	9.5%

样品类型 Cell culture supernatant, Serum, Plasma

检测类型 Sandwich (quantitative)

灵敏度 < 0.75 pg/ml

范围 0.78 pg/ml - 25 pg/ml

回收率 89 %

检测时间 4h 00m

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Human IL-12 ELISA kit High Sensitivity is designed for the quantitative measurement of IL-12 in Human serum, plasma, buffered solutions and cell culture medium.

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

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

平台 Microplate

性能

存放说明 Store at +4°C. Please refer to protocols.

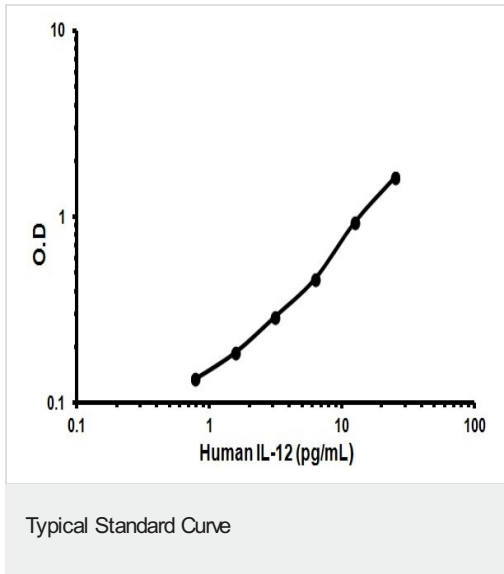
组件	标识符	1 x 96 tests	1 x 96 tests
10X Standard Diluent Buffer	Black	1 x 15ml	1 x 15ml
200X Wash Buffer	White	1 x 10ml	1 x 10ml
Amplification Diluent	Brown & blue spot	1 x 15ml	1 x 15ml
Amplifier	Yellow	1 x 200µl	1 x 200µl
Biotinylated Antibody Diluent	Red	1 x 7.5ml	1 x 7.5ml
Biotinylated anti-IL-12 p70	Red	1 x 0.4ml	1 x 0.4ml
Chromogen TMB Substrate Solution		1 x 11ml	1 x 11ml
HRP Diluent	Red	1 x 23ml	1 x 23ml
IL-12 Microplate (12 x 8 well strips)		1 unit	1 unit
IL-12 Standard (Lyophilized)	Yellow	2 vials	2 vials
Standard Diluent (Serum)		1 x 7ml	1 x 7ml
Stop Reagent	Black	1 x 11ml	1 x 11ml
Streptavidin-HRP		2 x 5µl	2 x 5µl

功能 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.

序列相似性 Belongs to the IL-6 superfamily.

细胞定位 Secreted.

图片



Representative Standard Curve using ab46035

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.cn/abpromise> or contact our technical team.

Terms and conditions

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