

# Human IL-12 p70 ELISA Kit ab100552

## 2 图像

### 概述

产品名称	人IL-12 p70 ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 1 pg/ml
范围	0.8 pg/ml - 600 pg/ml
回收率	88 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	95.52	88% - 109%
Plasma	85.63	79% - 101%

实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	Abcam's IL-12 p70 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human IL-12 p70 in plasma and cell culture supernatants.

This assay employs an antibody specific for Human IL-12 p70 coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-12 p70 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human IL-12 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-12 p70 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

We have not been able to detect endogenous Human IL-12 p70 in normal serum with ab100552, only in serum spiked with Human IL-12 p70.

Get results in 90 minutes with Human IL-12p40/p70 ELISA Kit ([ab178009](#)) from our SimpleStep ELISA® range.

**说明**

Optimization may be required with urine samples.  
The standard is affinity purified.

**平台**

Microplate

**性能**

**存放说明**

Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
300X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human IL-12 p70	2 vials
IL-12 p70 Microplate (12 x 8 wells)	1 unit
Recombinant IL-12 p70 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

**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.

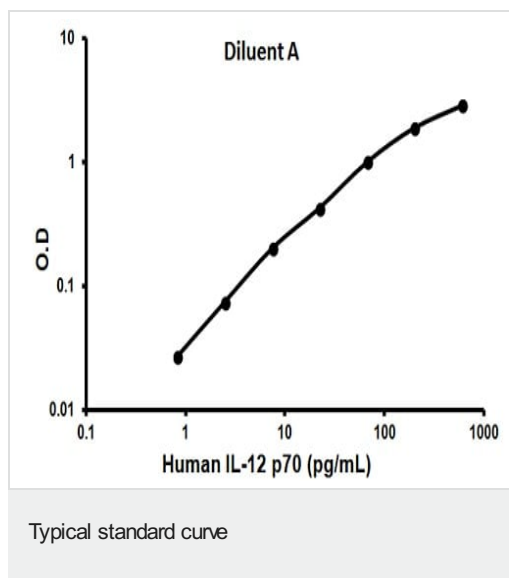
**序列相似性**

Belongs to the IL-6 superfamily.

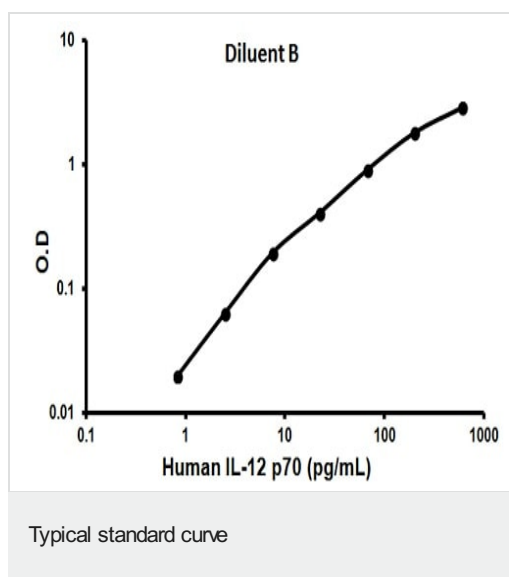
**细胞定位**

Secreted.

**图片**



Representative standard curve using ab100552



Representative standard curve using ab100552

**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