

Human IL-12 ELISA Kit High Sensitivity ab46035

5 References 画像数 1

製品の概要

製品名	Human IL-12 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
	Sample A	8	16.1pg/ml	1.25
	Sample B	8	6.1pg/ml	0.4
			CV%	
			7.8%	
			6.6%	
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Sample A	20	16pg/ml	1.2
	Sample B	20	6pg/ml	0.6
			CV%	
			7.9%	
			9.5%	
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	< 0.75 pg/ml			
検出範囲	0.78 pg/ml - 25 pg/ml			
添加回収試験	89 %			
全工程の試験時間	4h 00m			
ステップ	Multiple steps standard assay			
種交差性	交差種: Human			
製品の概要	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.

## 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	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

### 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.

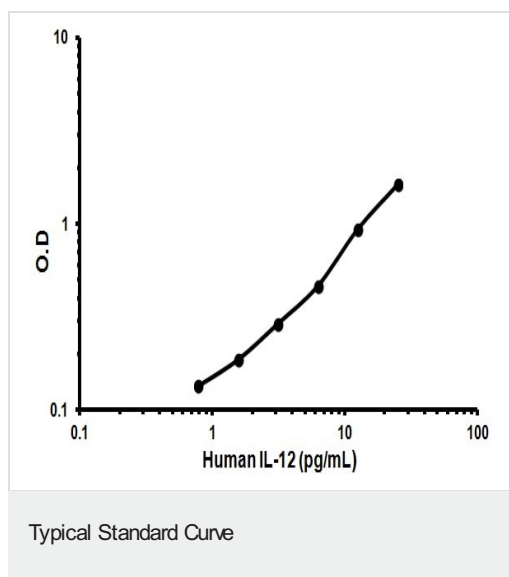
### 配列類似性

Belongs to the IL-6 superfamily.

### 細胞内局在

Secreted.

## 画像



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.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