

Human IL-12 p70 ELISA Kit ab46143

1 图像

概述				
产品名称	人IL-12 p70 ELISA试剂盒			
检测方法	Colorimetric			
精确度	批次内			
	样品	n	Mean	SD
	Overall			CV%
	批次间			
	样品	n	Mean	SD
	Overall	6		CV%
样品类型	Cell culture supernatant, Serum, Plasma			
检测类型	Sandwich (quantitative)			
灵敏度	= 2.2 pg/ml			
范围	6.25 pg/ml - 200 pg/ml			
回收率	特定样本回收率			
	样品类型	平均%		范围
	Serum	91		86% - 94%
检测时间	3h 40m			
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 (Interleukin- 12 p70) Human <i>in vitro</i> 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.

平台 Microplate

性能

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

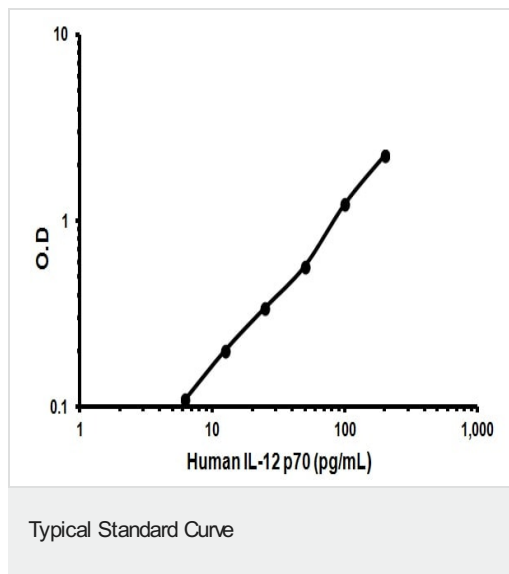
组 件	标识符	2 x 96 tests	1 x 96 tests
10X Standard Diluent Buffer	Black	1 x 25ml	1 x 15ml
200X Wash Buffer	White	2 x 10ml	1 x 10ml
Biotinylated Antibody Diluent	Red	1 x 13ml	1 x 7.5ml
Biotinylated anti-IL-12 p70	Red	2 x 400µl	1 x 400µl
Chromogen TMB Substrate Solution		1 x 24ml	1 x 11ml
Control	Silver	4 vials	2 vials
HRP Diluent	Red	1 x 23ml	1 x 12ml
IL-12 p70 Microplate (12 x 8 well strips)		2 units	1 unit
IL-12p70 Standard (Lyophilized)	Yellow	4 vials	2 vials
Standard Diluent (Serum)		2 x 7ml	1 x 7ml
Stop Reagent	Black	2 x 11ml	1 x 11ml
Streptavidin-HRP		4 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 ab46143

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