

Human IL-12 p70 ELISA Kit ab213791

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	CV%
1	16	18pg/ml	1.31	7.3%
2	16	51pg/ml	3.41	6.7%
3	16	294pg/ml	17.38	5.2%

批次间

样品	n	Mean	SD	CV%
1	24	19pg/ml	1.67	8.8%
2	24	55pg/ml	3.79	6.9%
3	24	310pg/ml	21.39	6.9%

样品类型 Cell culture supernatant, Serum

检测类型 Sandwich (quantitative)

灵敏度 < 2 pg/ml

范围 7.8 pg/ml - 500 pg/ml

检测时间 3h 30m

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 The Human IL-12(p70) Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213791) is designed for the quantitative detection of Human IL-12(p70) in cell culture supernatants and serum.

The ELISA kit is based on standard sandwich enzyme-linked immunosorbent assay technology. A

monoclonal antibody from mouse specific for IL-12(p70) has been pre-coated onto 96-well plates. Standards (Expression system for standard: sf21; Immunogen sequence: (I23-S328)+(R23-S219)) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for IL-12(p70) is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex was added and unbound conjugates were washed away with PBS or TBS buffer. HRP substrate TMB was used to visualize HRP enzymatic reaction. TMB was catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the Human IL-12(p70) amount of sample captured in plate. The standard product used in this kit is recombinant human IL-12(p70) with the molecular mass of 75 kDa.

说明

Interleukin (IL)-12 is a 70 kDa cytokine comprised of two disulfide-linked proteins (p35 and p40) and is essential for the initiation of effective immune response. The IL-12p70 is a heterodimer of p35 and p40 subunits; it is an important cytokine secreted by antigen-presenting cells in response to antigenic stimulation. Gene expression analysis of the IL-12 cytokine family subunits revealed that both strains induced high levels of p40 (protein chain communal to IL-12 p70 and IL-23) as well as p19, a subunit of IL-23. Conversely only ACT-18HS19 infection induced consistent transcription of IL-12 p35, a subunit of IL-12(p70).

平台

Pre-coated microplate (12 x 8 well strips)

性能

存放说明

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

组件	标识符	1 x 96 tests	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml	1 x 12ml
Adhesive Plate Seal		4 units	4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml	1 x 12ml
Anti-Human IL-12(p70) coated Microplate (12 x 8 wells)		1 unit	1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated anti- Human IL-12(p70) antibody		1 x 100µl	1 x 100µl
Lyophilized recombinant Human IL-12(p70) standard		2 vials	2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml	1 x 10ml
Wash Buffer (25X)		1 x 20ml	1 x 20ml

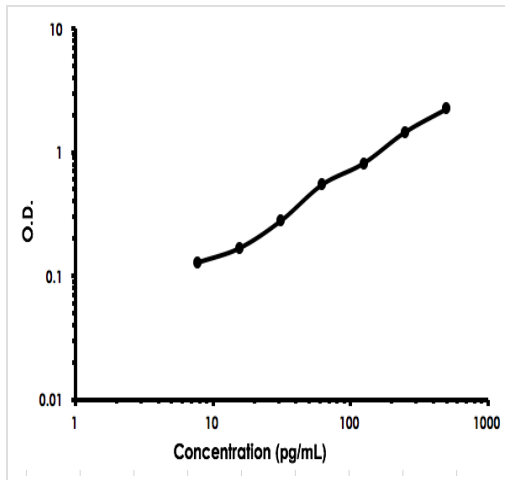
功能

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.

图片



Human IL-12(p70) ELISA Kit (ab213791) Standard Curve

Human IL-12(p70) ELISA Kit (ab213791) Standard
Curve

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