

Human IL-12 p70 ELISA Kit ab46143

[1 Image](#)

Overview

Product name Human IL-12 p70 ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall				4.3%

Inter-assay

Sample	n	Mean	SD	CV%
Overall	6			8.9%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity = 2.2 pg/ml

Range 6.25 pg/ml - 200 pg/ml

Recovery 95 %

Sample specific recovery

Sample type	Average %	Range
Serum	91	86% - 94%

Assay time 3h 45m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

Abcam's IL-12 p70 (Interleukin- 12 p70) Human *in vitro* 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.

Platform Microplate

Properties

Storage instructions Store at +4°C. Please refer to protocols.

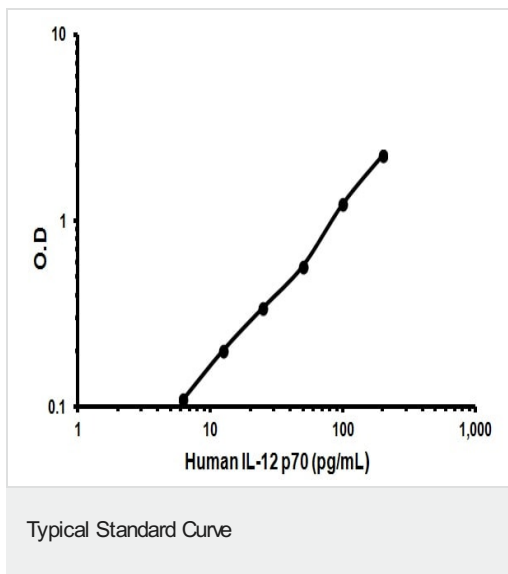
Components	Identifier	2 x 96 tests	1 x 96 tests
10X Standard Diluent Buffer	Black	1 x 25ml	1 x 15ml
200X Wash Buffer	White	2 x 10ml	1 x 10ml
Biotinylated Antibody Diluent	Red	1 x 13ml	1 x 7.5ml
Biotinylated anti-IL-12 p70	Red	2 x 400µl	1 x 400µl
Chromogen TMB Substrate Solution		1 x 24ml	1 x 11ml
Control	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

Function 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.

Sequence similarities Belongs to the IL-6 superfamily.

Cellular localization Secreted.

Images



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.com/abpromise> or contact our technical team.

Terms and conditions

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