

Product datasheet

Human IL-12 p70 ELISA Kit (Interleukin-12 p70)  
ab119532

1 Image

Overview

**Product name** Human IL-12 p70 ELISA Kit (Interleukin-12 p70)

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
Overall	8			3%

Inter-assay

Sample	n	Mean	SD	CV%
Overall	8			4.8%

**Sample type** Cell culture supernatant, Serum, Plasma

**Assay type** Sandwich (quantitative)

**Sensitivity** 2.1 pg/ml

**Range** 3.1 pg/ml - 200 pg/ml

**Recovery** 103 %

Sample specific recovery

Sample type	Average %	Range
Serum	103	79% - 118%

**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 accurate quantitative measurement of Human IL-12 p70 concentrations in Cell culture supernatant serum, and plasma (EDTA, citrate, heparin).

Human IL-12 p70 specific antibodies have been precoated onto 96-well plates. Standards and

test samples are added to the wells along with a biotin-conjugated Human IL-12 p70 detection antibody then incubated at room temperature. Following washing, a Streptavidin-HRP conjugate is added to each well, incubated at room temperature and washed. TMB is added and then catalyzed by HRP to produce a blue color product that changes into yellow after addition of acidic stop solution. The density of yellow coloration is directly proportional to the amount of Human IL-12 p70 captured on the plate.

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

**Tested applications**

**Suitable for:** Sandwich ELISA

**Platform**

Microplate

**Properties**

**Storage instructions**

Please refer to protocols.

Components	1 x 96 tests
20X Assay Buffer Concentrate	1 x 5ml
20X Wash Buffer Concentrate	1 x 50ml
Adhesive Films	4 units
Biotin-Conjugate anti-human IL12 p70 monoclonal antibody	1 x 100µl
Control high	1 vial
Control low	1 vial
IL-12 p70 Standard, lyophilized (400 pg/mL on reconstitution)	2 vials
Microplate coated with antibody to Human IL-12 p70 (12 x 8 wells)	1 unit
Sample Diluent	1 x 12ml
Stop Solution (1M Phosphoric acid)	1 x 15ml
Streptavidin-HRP	1 x 150µl
TMB Substrate Solution	1 x 15ml

**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.

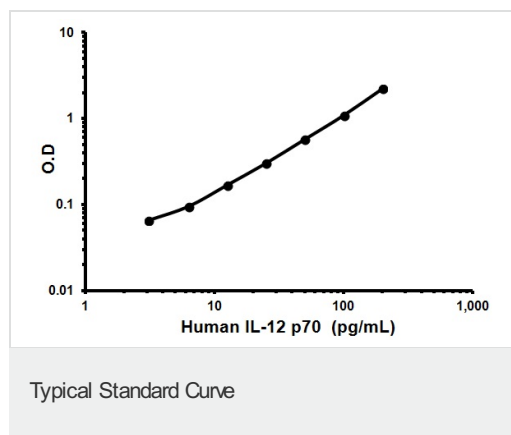
**Applications**

Our [Abpromise guarantee](#) covers the use of **ab119532** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

## Images



Representative Standard Curve using ab119532.

**Please note:** All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

## 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