

Mouse IL-12 p70 ELISA Kit ab119531

★★★★★ [1 Abreviews](#) [5 References](#) [1 Image](#)

Overview

Product name	Mouse IL-12 p70 ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall	7			8.3%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall	7			11%
Sample type	Cell culture supernatant, Serum				
Assay type	Sandwich (quantitative)				
Sensitivity	10 pg/ml				
Range	15.6 pg/ml - 1000 pg/ml				
Recovery	106 %				
	Sample specific recovery				
	Sample type		Average %		Range
	Serum		106		95% - 118%
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Mouse				
Product overview	IL-20 p70 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Mouse IL-20 p70 in serum and cell culture supernatants.				

IL-12 p70 specific antibodies have been precoated onto 96-well plates. Standards and test samples are added to the wells and along with a Biotin-conjugated anti-mouse IL-12 p70 detection antibody and the microplate is then incubated at room temperature. Following washing with wash buffer a Streptavidin-HRP conjugate is added to each well, incubated at room

temperature and then unbound conjugates are washed away using wash buffer. TMB is added and then catalyzed by HRP to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of IL-12 p70 captured on the plate.

**Notes** The standard is affinity purified.

**Platform** Microplate

## Properties

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

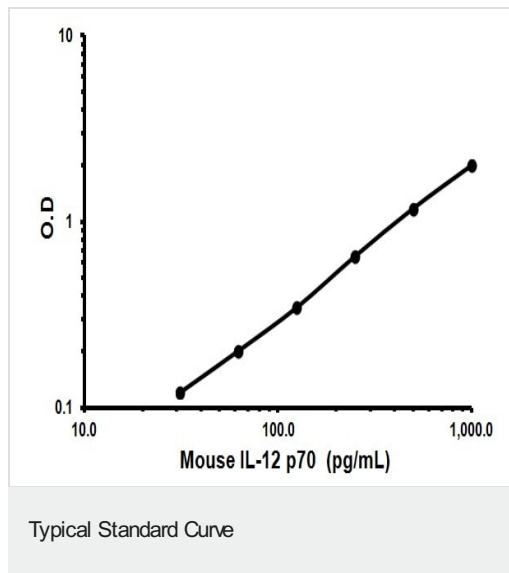
Components	1 x 96 tests
20X Assay Buffer Concentrate	1 x 5ml
20X Wash Buffer Concentrate	1 x 50ml
Adhesive Films	4 units
Biotin-Conjugate anti-mouse IL-12 p70 monoclonal antibody	1 x 70µl
Microplate coated with anti-mouse IL-12 p70 monoclonal antibody (12 x 8 wells)	1 unit
Mouse IL-12 p70 Standard, lyophilized (2 ng/mL on reconstitution).	2 vials
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.

## Images



Representative Standard Curve using ab119531.

**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