

Mouse IL-12 p70 ELISA Kit ab119531

★★★★★ [1 Abreviews](#) [5 References](#) [画像数 1](#)

製品の概要

製品名	Mouse IL-12 p70 ELISA Kit			
検出方法	Colorimetric			
再現性	Intra-Assay (同時再現性)			
	サンプル	N	平均値	SD
	Overall	7		CV%
				8.3%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Overall	7		CV%
				11%
サンプルの種類	Cell culture supernatant, Serum			
アッセイタイプ	Sandwich (quantitative)			
検出感度	10 pg/ml			
検出範囲	15.6 pg/ml - 1000 pg/ml			
添加回収試験	106 %			

	特定サンプルでの回収試験	
サンプルの種類	平均 %	測定範囲
Serum	106	95% - 118%

ステップ	Multiple steps standard assay
種交差性	交差種: Mouse
製品の概要	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.

**特記事項** The standard is affinity purified.

**試験プラットフォーム** Microplate

## 製品の特性

**保存方法** Store at +4°C. Please refer to protocols.

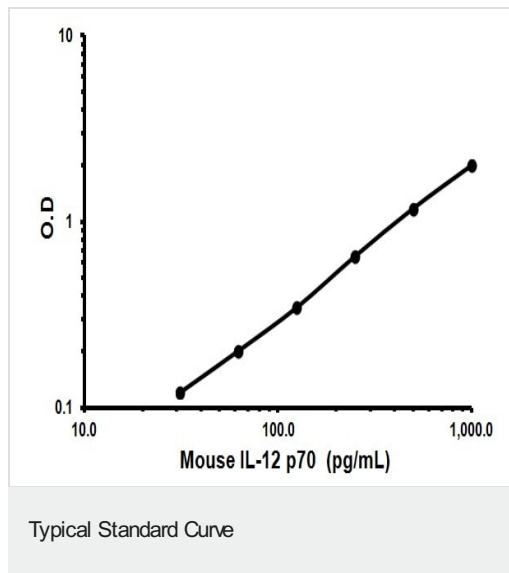
内容	1 x 96 tests
20X Assay Buffer Concentrate	1 x 5ml
20X Wash Buffer Concentrate	1 x 50ml
Adhesive Films	4 units
Biotin-Conjugate anti-mouse IL-12 p70 monoclonal antibody	1 x 70µl
Microplate coated with anti-mouse IL-12 p70 monoclonal antibody (12 x 8 wells)	1 unit
Mouse IL-12 p70 Standard, lyophilized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

**機能** 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.

**配列類似性** Belongs to the IL-6 superfamily.

**細胞内局在** Secreted.

## 画像



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.co.jp/abpromise> or contact our technical team.

### Terms and conditions

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