

Monkey IL-2 ELISA Kit ab119541

画像数 1

製品の概要

製品名	Monkey IL-2 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	8		CV%
				< 7%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Overall	8		CV%
				< 5%
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 U/ml			
検出範囲	20 U/ml - 1250 U/ml			
添加回収試験	80 %			
ステップ	Multiple steps standard assay			
種交差性	交差種: Monkey			
製品の概要	Abcam's IL-2 Monkey <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for accurate quantitative measurement of Monkey IL-2 concentrations in cell culture supernatant, serum and plasma (EDTA, citrate, heparin), amniotic fluid, urine and whole blood.			

IL-2 specific antibodies have been precoated onto 96-well plates. Standards and test samples are added to the wells along with a biotin-conjugated IL-2 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 the addition of an acidic stop solution. The density of yellow coloration is directly proportional to the amount of IL-2 captured on the plate.

特記事項 Standard is recombinant protein produced in *E.coli*.

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-monkey IL-2 polyclonal antibody	1 x 70µl
Microplate coated with monoclonal antibody to monkey IL-2	1 unit
Monkey IL-2 Standard lyophilized	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

Produced by T-cells in response to antigenic or mitogenic stimulation, this protein is required for T-cell proliferation and other activities crucial to regulation of the immune response. Can stimulate B-cells, monocytes, lymphokine-activated killer cells, natural killer cells, and glioma cells.

関連疾患

Note=A chromosomal aberration involving IL2 is found in a form of T-cell acute lymphoblastic leukemia (T-ALL). Translocation t(4;16)(q26;p13) with involves TNFRSF17.

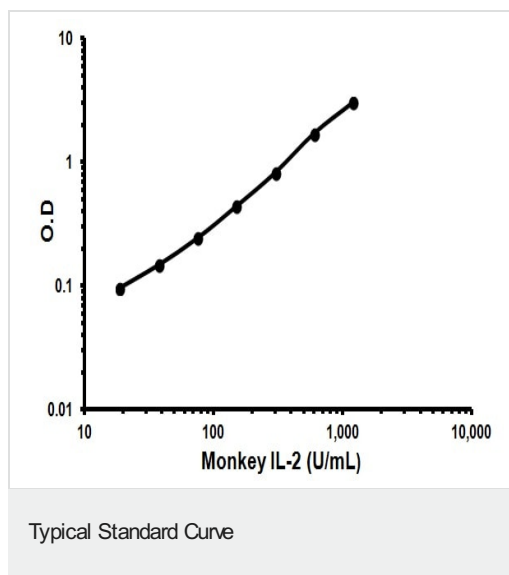
配列類似性

Belongs to the IL-2 family.

細胞内局在

Secreted.

画像



Representative Standard Curve using ab119541.

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