



Product datasheet

Rat IL-17A ELISA Kit ab119536

4 References 画像数 1

製品の概要

製品名	Rat IL-17A 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%
				8.5%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Overall	7		CV%
				7.6%
サンプルの種類	Cell culture supernatant, Serum, Plasma			
アッセイタイプ	Sandwich (quantitative)			
検出感度	1 pg/ml			
検出範囲	1.6 pg/ml - 100 pg/ml			
添加回収試験	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
	Cell culture supernatant			92% - 112%
	Serum			34% - 44%

ステップ	Multiple steps standard assay
種交差性	交差種: Rat
製品の概要	Abcam's IL-17A Rat <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for accurate quantitative measurement of Rat IL-17 (Interleukin-17) concentrations in Cell culture supernatant, serum and plasma (EDTA, citrate, heparin).

IL-17A specific antibodies have been precoated onto 96-well plates. Standards and test samples are added to the wells along with a biotin-conjugated IL-17A 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 IL-17A captured on the plate.

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-rat IL-17A monoclonal antibody	1 x 70µl
Microplate coated with monoclonal antibody to rat IL-17A (12 x 8 wells)	1 unit
Rat IL-17A Standard Lyophilized (200 pg/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

Induces stromal cells to produce proinflammatory and hematopoietic cytokines. Enhances the surface expression of the intracellular adhesion molecule-1 (ICAM-1) in fibroblasts.

組織特異性

Restricted to activated memory T-cells.

配列類似性

Belongs to the IL-17 family.

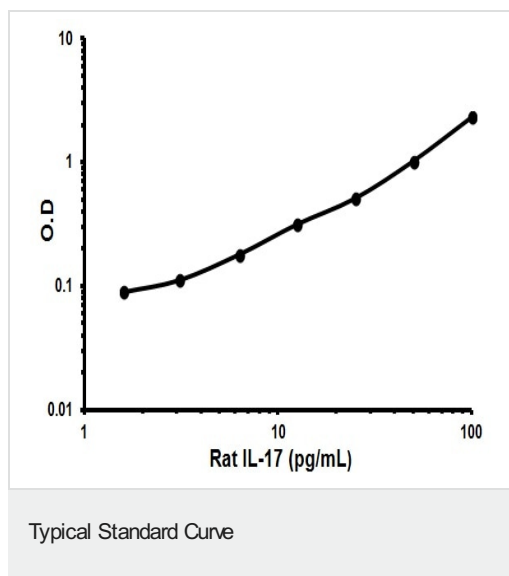
翻訳後修飾

Found both in glycosylated and nonglycosylated forms.

細胞内局在

Secreted.

画像



Representative Standard Curve using ab119536.

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