

Rat IL-17F ELISA Kit ab213913

画像数 1

製品の概要

製品名	Rat IL-17F 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	CV%
	Sample 1	16	275pg/ml	12.93	4.7%
	Sample 2	16	524pg/ml	32.5	6.2%
	Sample 3	16	928pg/ml	52	5.6%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	Sample 1	24	352pg/ml	17.6	5%
	Sample 2	24	723pg/ml	52.78	7.3%
	Sample 3	24	1230pg/ml	83.64	6.8%
サンプルの種類	Cell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma, Cit plasma, Tissue Homogenate				
アッセイタイプ	Sandwich (quantitative)				
検出感度	< 10 pg/ml				
検出範囲	31.2 pg/ml - 2000 pg/ml				
全工程の試験時間	3h 30m				
ステップ	Multiple steps standard assay				
種交差性	交差種: Rat				
製品の概要	The Rat IL-17F Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213913) is designed for the quantitative measurement of Rat IL-17F in cell culture supernatants, cell lysates, tissue homogenates, serum and plasma (heparin, EDTA).				

The ELISA kit is based on standard sandwich enzyme-linked immune-sorbent assay technology. A monoclonal antibody from mouse specific for IL-17F has been pre-coated onto 96-well plates. Standards and test samples are added to the wells; a biotinylated detection polyclonal antibody from goat specific for IL-17F is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. HRP substrate TMB is used to visualize HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the rat IL-17F amount of sample captured in plate.

特記事項

Interleukin 17F, also called IL17F is involved in the regulation of normal versus aberrant T-cell responses. This gene is mapped to 6p12.2. The protein encoded by this gene is a cytokine that shares sequence similarity with IL17. This cytokine is expressed by activated T cells, and has been shown to stimulate the production of several other cytokines, including IL6, IL8, and CSF2/GM-CSF. This cytokine is also found to inhibit the angiogenesis of endothelial cells and induce endothelial cells to produce IL2, TGFB1/TGFB, and monocyte chemoattractant protein-1. It is suggested that targeting IL17 and IL17F or antagonizing IL17R might mitigate neutrophil-mediated inflammation in CF.

試験プラットフォーム

Pre-coated microplate (12 x 8 well strips)

製品の特性

保存方法

Store at -20°C. Please refer to protocols.

内容	ラベル	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Adhesive Plate Seal		4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-rat IL-17F coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti- rat IL-17F antibody		1 x 100µl
Lyophilized recombinant rat IL-17F standard		2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml
Wash Buffer (25X)		1 x 20ml

機能

Ligand for IL17RA and IL17RC (PubMed:17911633). The heterodimer formed by IL17A and IL17F is a ligand for the heterodimeric complex formed by IL17RA and IL17RC (PubMed:18684971). Involved in stimulating the production of other cytokines such as IL6, IL8 and CSF2, and in regulation of cartilage matrix turnover (PubMed:11591732, PubMed:11591768,

PubMed:11574464). Also involved in stimulating the proliferation of peripheral blood mononuclear cells and T-cells and in inhibition of angiogenesis (PubMed:11591732). Plays a role in the induction of neutrophilia in the lungs and in the exacerbation of antigen-induced pulmonary allergic inflammation.

組織特異性

Expressed in activated, but not resting, CD4+ T-cells and activated monocytes.

関連疾患

Candidiasis, familial, 6

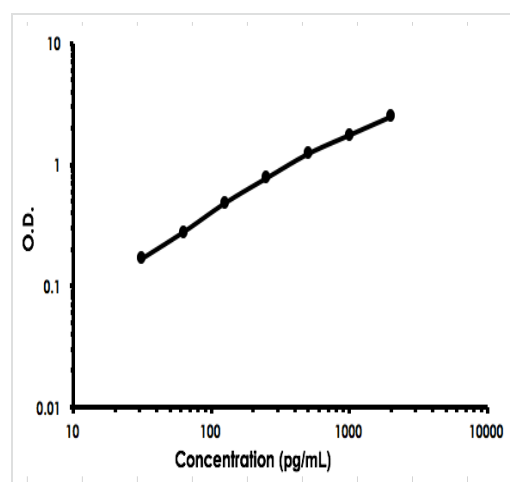
配列類似性

Belongs to the IL-17 family.

細胞内局在

Secreted.

画像



Rat IL-17F ELISA Kit (ab213913) Standard Curve

Rat IL-17F ELISA Kit (ab213913) Standard Curve

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