

Rat IL-18 ELISA Kit ab213909

14 References 画像数 1

製品の概要

製品名	Rat IL-18 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	29pg/ml	1.97	6.8%
	Sample 2	16	166pg/ml	2.2	4.7%
	Sample 3	16	511pg/ml	22.99	4.5%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	Sample 1	24	29pg/ml	2.2	7.6%
	Sample 2	24	165pg/ml	9.07	5.5%
	Sample 3	24	490pg/ml	29.4	6%
サンプルの種類	Cell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma				
アッセイタイプ	Sandwich (quantitative)				
検出感度	< 1 pg/ml				
検出範囲	15.6 pg/ml - 1000 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-18 Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213909) is designed for the				

The ELISA kit is based on standard sandwich enzyme-linked immune-sorbent assay technology. A monoclonal antibody from mouse specific for IL-18 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-18 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-18 amount of sample captured in plate.

特記事項

Interleukin (IL)-18, also called Interferon-gamma-inducing factor (IGIF), augments natural killer(NK) activity in spleen cells. The gene encodes a precursor protein of 192 amino acids and a mature protein of 157 amino acids. IL-18 is a recently discovered cytokine that modulates both T helper type 1 (Th1) and Th2 responses. IL-18 is a potent proinflammatory cytokine with potential atherogenic properties. It is highly expressed in the atherosclerotic plaques compared with control normal arteries and is localized mainly in plaque macrophages.

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	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml	1 x 12ml
Adhesive Plate Seal		4 units	4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml	1 x 12ml
Anti-rat IL-18 coated Microplate (12 x 8 wells)		1 unit	1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated anti- Rat IL-18 antibody		1 x 100µl	1 x 100µl
Lyophilized recombinant Rat IL-18 standard		2 vials	2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml	1 x 10ml
Wash Buffer (25X)		1 x 20ml	1 x 20ml

機能

Augments natural killer cell activity in spleen cells and stimulates interferon gamma production in T-helper type I cells.

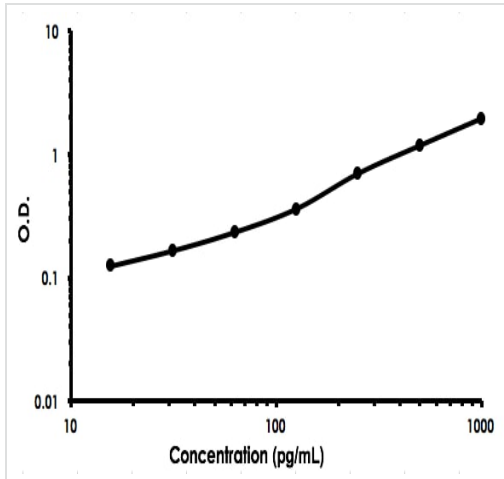
配列類似性

Belongs to the IL-1 family.

細胞内局在

Secreted.

画像



Rat IL-18 ELISA Kit (ab213909) Standard Curve

Rat IL-18 ELISA Kit (ab213909) 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