

Rat IL-10 ELISA Kit ab108872

3 References 画像数 1

製品の概要

製品名	Rat IL-10 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			CV%
				4.3%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				9.9%
サンプルの種類	Cell culture supernatant			
アッセイタイプ	Sandwich (quantitative)			
検出感度	= 45 pg/ml			
検出範囲	0.063 ng/ml - 4 ng/ml			
全工程の試験時間	4h 30m			
ステップ	Multiple steps standard assay			
種交差性	交差種: Rat			
製品の概要	Abcam's IL-10 Rat <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-10 in cell culture supernatants.			

An IL-10 specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently an IL-10 specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase 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-10 captured in plate.

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

製品の特性**保存方法**

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

内容	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
1X Standard Diluent	1 x 2ml
20X Wash Buffer Concentrate	2 x 30ml
60X Biotinylated Rat IL-10 Antibody	1 x 100µl
Chromogen Substrate	1 x 7ml
IL-10 Microplate (12 x 8 well strips)	1 unit
IL-10 Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

機能

Inhibits the synthesis of a number of cytokines, including IFN-gamma, IL-2, IL-3, TNF and GM-CSF produced by activated macrophages and by helper T-cells.

組織特異性

Produced by a variety of cell lines, including T-cells, macrophages, mast cells and other cell types.

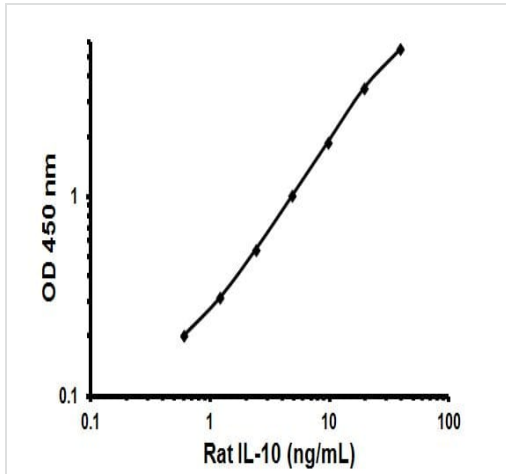
配列類似性

Belongs to the IL-10 family.

細胞内局在

Secreted.

画像



Representative Standard Curve Using ab108872.

Typical 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