

Rat IL-10 ELISA Kit ab214566

リコンビナント SimpleStep ELISA

21 References 画像数 7

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	CV%
	Overall	8			5.6%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	Overall	3			3.4%
サンプルの種類	Cell culture supernatant, Serum, Cit plasma				
アッセイタイプ	Sandwich (quantitative)				
検出感度	4.89 pg/ml				
検出範囲	31.3 pg/ml - 2000 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	93		84% - 109%	
	Serum	94		93% - 94%	
	Cit plasma	86		85% - 89%	
全工程の試験時間	1h 30m				
ステップ	One step assay				
種交差性	交差種: Rat 非交差種: Cow				
製品の概要	Rat IL-10 ELISA Kit (ab214566) is a single-wash 90 min sandwich ELISA designed for the				

quantitative measurement of IL-10 protein in cell culture supernatant, cit plasma, and serum. It uses our proprietary SimpleStep ELISA® technology. Quantitate Rat IL-10 with 4.89 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate (**ab203359**) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

特記事項

Interleukin-10 is an anti-inflammatory cytokine involved in inhibition of a variety of other cytokines, including IFN-gamma, TNF, IL-2 and GM-CSF.

試験プラットフォーム

Pre-coated microplate (12 x 8 well strips)

製品の特性

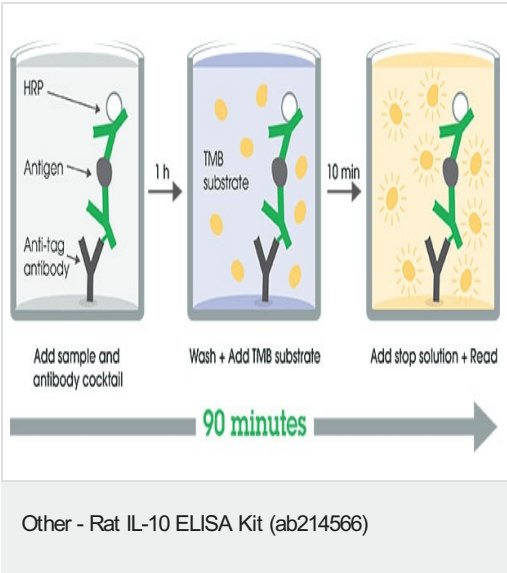
保存方法

Store at +4°C. Please refer to protocols.

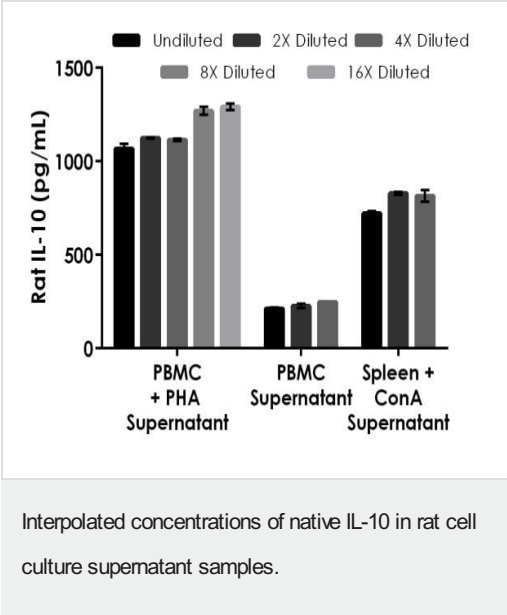
内容	1 x 96 tests
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 5BR	1 x 6ml
Plate Seals	1 unit
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
10X Rat IL-10 Capture Antibody	1 x 600µl
10X Rat IL-10 Detector Antibody	1 x 600µl
Rat IL-10 Lyophilized Recombinant Protein	2 vials
Sample Diluent 25BS	1 x 20ml
Sample Diluent NS (ab193972)	1 x 50ml
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

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.

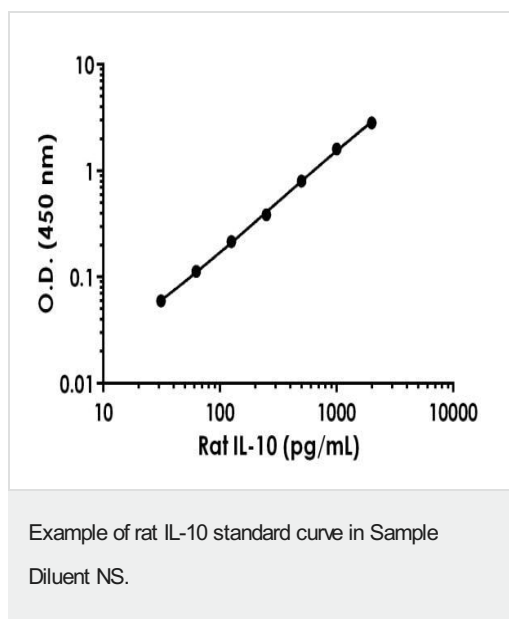
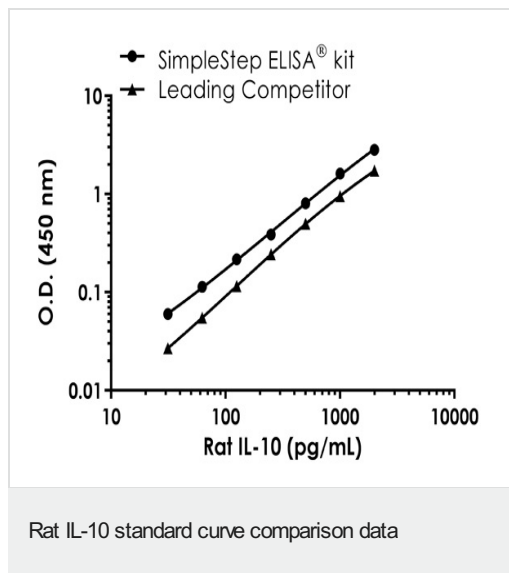
画像

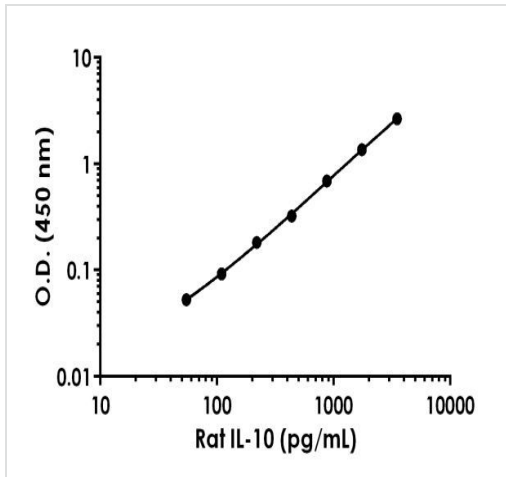


SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



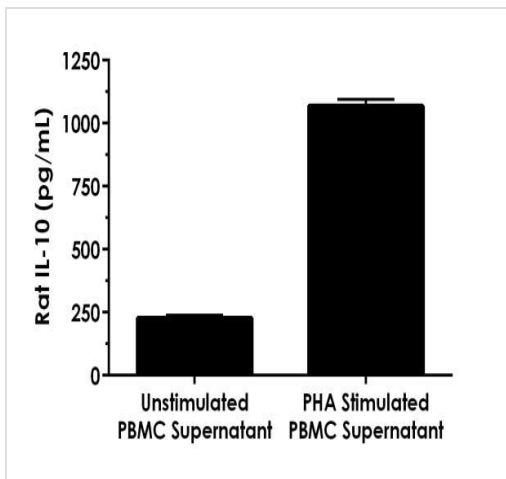
The concentrations of IL-10 were measured in duplicates, interpolated from the IL-10 standard curves and corrected for sample dilution. Undiluted samples are as follows: rat PBMC stimulated with PHA (10 µg/mL, 48 hours) supernatant 50%, unstimulated PBMC supernatant 100%, and spleen stimulated with concanavalin A (5 µg/mL, 48 hours) supernatant 50%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean IL-10 concentration was determined to be 1,174 pg/mL in PBMC stimulated with PHA supernatant, 230 pg/mL in unstimulated PBMC supernatant and 789 pg/mL in spleen stimulated with concanavalin A supernatant.





Example of rat IL-10 standard curve in Sample Diluent 25BS.

Background-subtracted data values (mean \pm SD) are graphed.



Rat peripheral blood mononuclear cells were cultured unstimulated or stimulated with 10 μ g/mL PHA.

Conditioned media was harvested after 48 hours. IL-10 was measured in 50% unstimulated and PHA stimulated PBMC supernatant. The concentrations of IL-10 were measured in duplicate, interpolated from the IL-1beta standard curves and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean \pm SD, n=2). The mean IL-10 concentration was determined to be 1,068 pg/mL in PHA stimulated PBMC supernatant and 227 pg/mL in unstimulated PBMC supernatant.

Powered by
recombinant antibodies




Research with confidence
Consistent and reproducible results



Long-term and scalable supply
Recombinant technology



Success from the first experiment
Confirmed specificity



Ethical standards compliant
Animal-free production

Sandwich ELISA - Rat IL-10 ELISA Kit (ab214566)

To learn more about the advantages of recombinant antibodies see [here](#).

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