

Human IL-10 ELISA Kit High Sensitivity ab46059

5 References 画像数 1

製品の概要

製品名	Human IL-10 ELISA Kit High Sensitivity				
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	3			7.8%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	Overall	3			10.2%
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	< 0.98 pg/ml				
検出範囲	1.56 pg/ml - 50 pg/ml				
添加回収試験	87 %				
	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		94	% - %	
	Serum		80	59% - 93%	
全工程の試験時間	3h 00m				
ステップ	Multiple steps standard assay				
種交差性	交差種: Human				
製品の概要	Abcam's IL-10 Human High Sensitivity <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-10 in Human sera, plasmas, buffered solutions and cell culture media.				

A monoclonal antibody specific for IL-10 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IL-10 concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for IL-10 are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated and washed. A TMB substrate solution is added which acts on the bound enzyme to induce a colored reaction product. The intensity of this colored product is directly proportional to the concentration of IL-10 present in the samples.

This kit will recognize both endogenous and recombinant Human IL-10.

Get results in 90 minutes with Human IL-10 ELISA Kit ([**ab185986**](#)) from our SimpleStep ELISA® range.

試験プラットフォーム

Microplate

製品の特長

保存方法

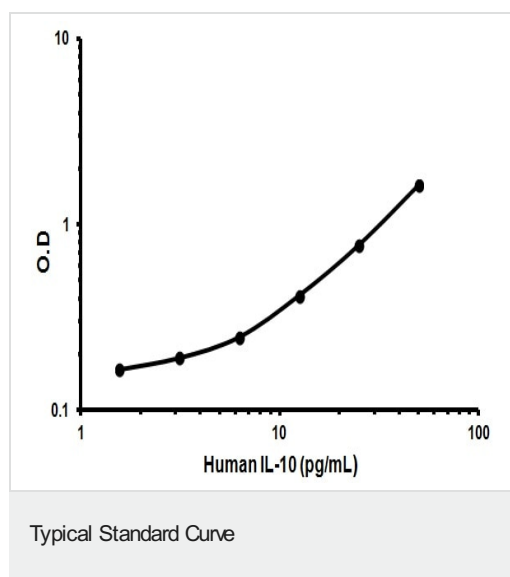
Store at +4°C. Please refer to protocols.

内容	ラベル	1 x 96 tests
10X Standard Diluent Buffer	Black	1 x 15ml
200X Wash Buffer	White	1 x 10ml
Amplification Diluent	Brown & blue spot	1 x 15ml
Amplifier	Yellow	1 x 200µl
Biotinylated anti IL-10	Red	1 x 400µl
Biotinylated Antibody Diluent	Red	1 x 7.5ml
Chromogen TMB Substrate Solution		1 x 11ml
HRP Diluent	Red	1 x 23ml
IL-10 Microplate (12 x 8 well strips)		1 unit
IL-10 Standard (Lyophilized)	Yellow	2 vials
Stop Reagent	Black	1 x 11ml

内容	ラベル	1 x 96 tests
Streptavidin-HRP		2 x 5µl

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 ab46059

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