

Human IL-6 Receptor ELISA Kit ab46029

★★★★★ 1 Abreviews 6 References 画像数 1

製品の概要

製品名	Human IL-6 Receptor 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 A	7	977.7pg/ml	3.04	0.31%
	Sample B	7	131.1pg/ml	2.489	2.289%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	Sample A	24	983.9pg/ml	10.46	1.06%
	Sample B	24	136.2pg/ml	6.086	4.46%
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	< 5 pg/ml				
検出範囲	31.25 pg/ml - 1000 pg/ml				
添加回収試験	96 %				
全工程の試験時間	1h 45m				
ステップ	Multiple steps standard assay				
種交差性	交差種: Human				
製品の概要	Abcam's IL-6 Receptor (Interleukin-6 Receptor) Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of soluble IL-6 Receptor in Human serum, plasma, buffered solutions and cell culture medium.				

A monoclonal antibody specific for IL-6 Receptor has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IL-6 Receptor 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-6 Receptor 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-6 Receptor present in the samples.

This kit will recognize both endogenous and recombinant Human IL-6 Receptor.

Get higher sensitivity in only 90 minutes with Human IL-6 Receptor ELISA Kit (**ab181424**) 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	2 x 96 tests
10X Standard Diluent Buffer	Black	1 x 15ml	1 x 25ml
200X Wash Buffer	White	1 x 10ml	2 x 10ml
Biotinylated Antibody Diluent	Red	1 x 7ml	1 x 13ml
Biotinylated anti-IL-6R (CD126)	Red	1 x 0.4ml	2 x 0.4ml
Chromogen TMB Substrate Solution		1 x 11ml	1 x 24ml
HRP Diluent	Red	1 x 12ml	1 x 23ml
IL-6 Receptor Microplate (12 x 8 well strips)		1 unit	2 units
IL-6R Control	Silver	2 vials	4 vials
IL-6R Standard	Yellow	2 vials	4 vials
Stop Reagent	Black	1 x 11ml	2 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl

機能

Part of the receptor for interleukin 6. Binds to IL6 with low affinity, but does not transduce a signal. Signal activation necessitate an association with IL6ST. Activation may lead to the regulation of the immune response, acute-phase reactions and hematopoiesis.
Low concentration of a soluble form of IL6 receptor acts as an agonist of IL6 activity.

組織特異性

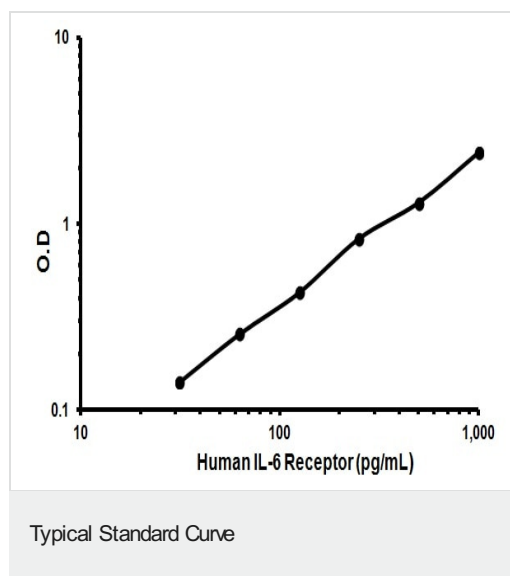
Isoform 2 is expressed in peripheral blood mononuclear cells and weakly found in urine and serum.

配列類似性

Belongs to the type I cytokine receptor family. Type 3 subfamily.
Contains 1 fibronectin type-III domain.
Contains 1 Ig-like C2-type (immunoglobulin-like) domain.

ドメイン	<p>The two fibronectin type-III-like domains, contained in the N-terminal part, form together a cytokine-binding domain.</p> <p>The WSXWS motif appears to be necessary for proper protein folding and thereby efficient intracellular transport and cell-surface receptor binding.</p>
翻訳後修飾	A short soluble form may also be released from the membrane by proteolysis.
細胞内局在	Secreted and Basolateral cell membrane.

画像



Representative Standard Curve using ab46029

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