

Human CD130 ELISA Kit (GP130) ab46135

1 References 画像数 1

製品の概要

製品名 Human CD130 ELISA Kit (GP130)

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%
A	8	1751pg/ml	17.61	1%
B	8	264.3pg/ml	8.67	3.27%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
A	16	1782pg/ml	26.64	1.49%
B	16	247pg/ml	11.99	4.85%

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 < 15 pg/ml

検出範囲 56.25 pg/ml - 1800 pg/ml

添加回収試験 97 %

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
Serum	97	% - %

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 CD130 (GP130) Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of CD130 in Human serum, plasma, buffered solutions and supernatants.

A monoclonal antibody specific for CD130 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known CD130 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 CD130 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 CD130 present in the samples.

This kit will recognize both endogenous and recombinant Human CD130.

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.

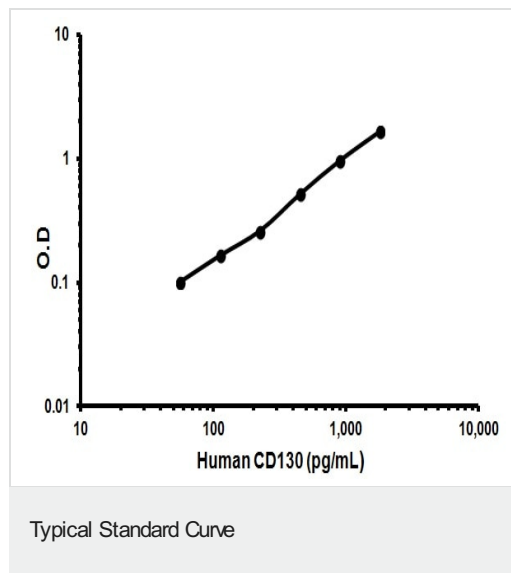
内容	ラベル	2 x 96 tests	1 x 96 tests	1 x 96 tests
10X Standard Diluent Buffer	Black	1 x 25ml	1 x 15ml	1 x 15ml
200X Wash Buffer	White	2 x 10ml	1 x 10ml	1 x 10ml
Biotinylated Antibody Diluent	Red	1 x 13ml	1 x 7.5ml	1 x 7.5ml
Biotinylated anti-CD130	Red	2 x 400µl	1 x 400µl	1 x 400µl
CD130 Microplate (12 x 8 well strips)		2 units	1 unit	1 unit
CD130 Standard (lyophilized)	Yellow	4 vials	2 vials	2 vials
Chromogen TMB Substrate Solution		1 x 24ml	1 x 11ml	1 x 11ml
Control	Silver	4 vials	2 vials	2 vials
HRP Diluent	Red	1 x 23ml	1 x 12ml	1 x 12ml
Stop Reagent	Black	2 x 11ml	1 x 11ml	1 x 11ml
Streptavidin-HRP		4 x 5µl	2 x 5µl	2 x 5µl

関連性

CD130 is a signal transducer shared by many cytokines, including interleukin 6 (IL6), ciliary neurotrophic factor (CNTF), leukemia inhibitory factor (LIF), and oncostatin M (OSM). This protein functions as a part of the cytokine receptor complex. The activation of this protein is dependent upon the binding of cytokines to their receptors. vIL6, a protein related to IL6 and encoded by the Kaposi sarcoma-associated herpesvirus, can bypass the interleukin 6 receptor (IL6R) and directly activate this protein. Knockout studies in mice suggested a critical role of the gene encoding this protein in regulating myocyte apoptosis. Alternatively spliced transcript variants encoding distinct isoforms have been described.

細胞内局在

Cell membrane; single-pass type I membrane protein and secreted



Representative Standard Curve using ab46135

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