

Human IP-10 ELISA Kit (CXCL10) ab83700

8 References 画像数 1

製品の概要

製品名 Human IP-10 ELISA Kit (CXCL10)

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	6	60.1pg/ml	0.9	1.6%
B	6	32.4pg/ml	0.5	1.5%
C	6	19.6pg/ml	0.3	1.3%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
A	18	56.9pg/ml	2.7	4.7%
B	18	29.9pg/ml	2	6.5%
C	18	17.7pg/ml	1.5	8.5%

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 < 5.7 pg/ml

検出範囲 6.25 pg/ml - 200 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
Serum	120	118.7% - 122.4%

全工程の試験時間 3h 45m

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要

Abcam’s IP-10 (CXCL10) Human in vitro ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IP-10 in supernatants, buffered solutions, serum and plasma samples.

A monoclonal antibody specific for IP-10 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IP-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 IP-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 IP-10 present in the samples.

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

試験プラットフォーム

Microplate

製品の特性

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
10X Standard Diluent Buffer	Black	1 x 25ml	1 x 15ml
200X Wash Buffer	White	2 x 10ml	1 x 10ml
Biotinylated anti IP-10		2 x 400µl	1 x 400µl
Biotinylated Antibody Diluent	Red	1 x 13ml	1 x 7ml
Chromogen TMB Substrate Solution		1 x 24ml	1 x 11ml
HRP Diluent	Red	1 x 23ml	1 x 12ml
IP-10 Microplate (12 x 8 well strips)		2 units	1 unit
IP-10 Standard (Lyophilized)		4 vials	2 vials
Plastic Plate Covers		4 units	2 units
Stop Reagent	Black	2 x 11ml	1 x 11ml
Streptavidin-HRP		4 x 5µl	2 x 5µl

機能

Chemotactic for monocytes and T-lymphocytes. Binds to CXCR3.

配列類似性

Belongs to the intercrine alpha (chemokine CxC) family.

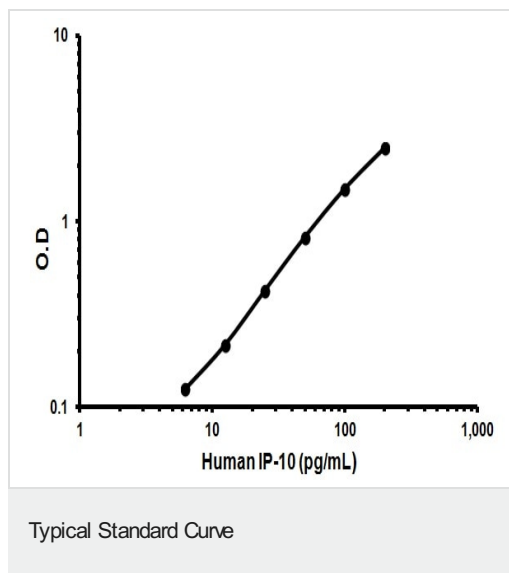
翻訳後修飾

CXCL10(1-73) is produced by proteolytic cleavage after secretion from keratinocytes.

細胞内局在

Secreted.

画像



Representative Standard Curve using ab83700

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