

Human CXCL16 ELISA Kit ab187397

SimpleStep ELISA

画像数 5

製品の概要

製品名	Human CXCL16 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%
Human serum	5			1.5%

サンプル	N	平均値	SD	CV%
Human serum	3			5.6%

サンプルの種類	Cell culture supernatant, Serum, Hep Plasma, EDTA Plasma, Cit plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	2 pg/ml
検出範囲	31.25 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
Serum	81.2	79.5% - 82.5%
Cell culture media	108	104.3% - 114.5%
Hep Plasma	87.1	73.5% - 101.3%
EDTA Plasma	87.9	80.8% - 91.8%
Cit plasma	87	85.9% - 87.8%

全工程の試験時間	1h 30m
ステップ	One step assay

## 種交差性

交差種: Human

## 製品の概要

Human CXCL16 ELISA Kit (ab187397) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of CXCL16 protein in cell culture supernatant, cit plasma, edta plasma, hep plasma, and serum. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human CXCL16 with 2 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

CXCL16 is a member of the chemokine superfamily. CXCL16 may act as a scavenger receptor on macrophages, which specifically binds to OxLDL (oxidized low density lipoprotein), suggesting that it may be involved in pathophysiology such as atherogenesis. CXCL16 induces a strong chemotactic response. CXCL16 may have a role in leukocytes migration. CXCL16 induces calcium mobilization.

## 試験プラットフォーム

Microplate

## 製品の特性

### 保存方法

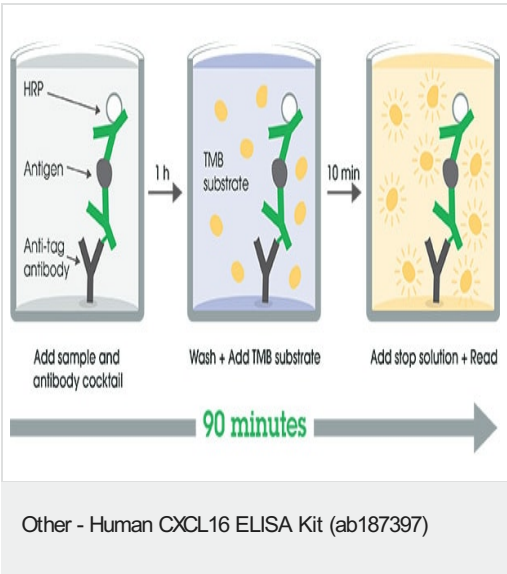
Store at +4°C. Please refer to protocols.

内容	1 x 96 tests
10X Human CXCL16 Capture Antibody	1 x 600µl
10X Human CXCL16 Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent CPI - HAMA Blocker (ab193969)	1 x 6ml
Human CXCL16 Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit

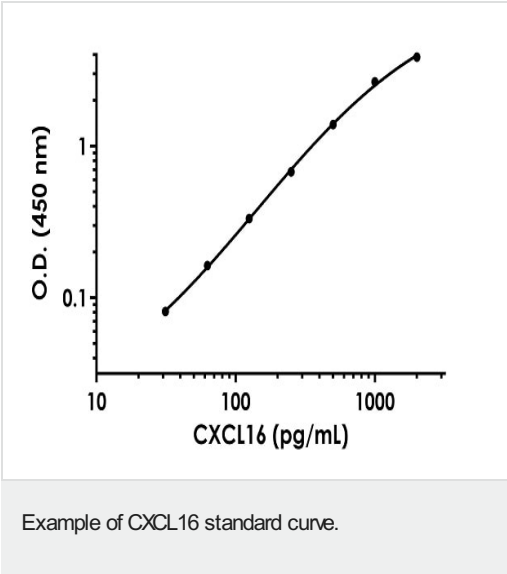
内容	1 x 96 tests
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

<div> <div>関連性</div> <div> <p>CXCL16 acts as a scavenger receptor on macrophages, which specifically binds to OxLDL (oxidized low density lipoprotein), suggesting that it may be involved in pathophysiology such as atherogenesis. It induces a strong chemotactic response and calcium mobilization. It binds to CXCR6/Bonzo.</p> </div> </div> <div> <div>細胞内局在</div> <div> <p>Cell Membrane and Secreted</p> </div> </div>
---

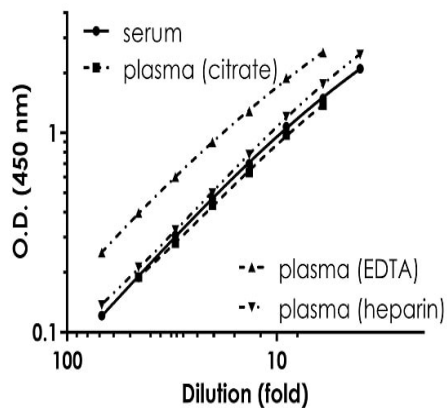
画像



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.

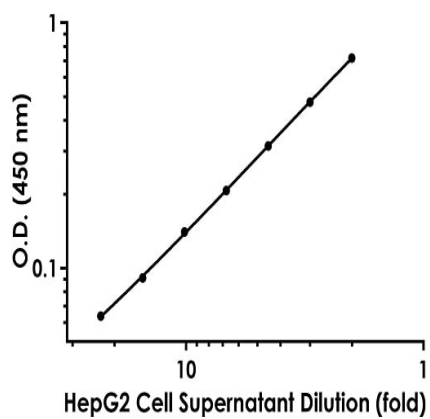


Background-subtracted data values (mean +/- SD) are graphed.



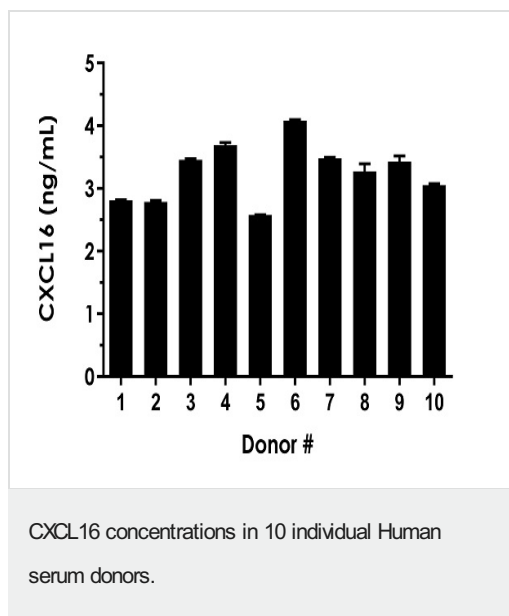
Background-subtracted data values (mean  $\pm$  SD,  $n = 2$ ) are graphed.

Titration of Human serum and plasma within the working range of the assay.



Background-subtracted data values (mean  $\pm$  SD,  $n = 2$ ) are graphed.

Titration of HepG2 cell culture supernatant within the working range of the assay.



8 X diluted serum samples from 10 apparently healthy male donors were measured using this kit. Interpolated data values corrected for sample dilution are graphed in ng of CXCL16 per mL of serum (mean  $\pm$  SD, n=3). The mean of CXCL16 concentration of these ten individual serum samples was determined to be 3.26 ng /mL with a range of 2.58 – 4.08 ng /mL.

**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