

### Human CXCL9 ELISA Kit, Fluorescent ab270891

リコンビナント SimpleStep ELISA

画像数 6

#### 製品の概要

製品名 Human CXCL9 ELISA Kit, Fluorescent

検出方法 Fluorescent

再現性 Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
Supernatant	3			3.4%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Supernatant	5			4.9%

サンプルの種類 Cell culture supernatant, Tissue Extracts

アッセイタイプ Sandwich (quantitative)

検出感度 2.6 pg/ml

検出範囲 2.93 pg/ml - 3000 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
Cell culture supernatant	90	86% - 92%
Tissue Extracts	109	104% - 116%

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 CXCL9 *in vitro* CatchPoint® SimpleStep ELISA® (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of CXCL9 protein in human plasma cell culture supernatants, and tissue extract samples.

This CatchPoint SimpleStep ELISA kit has been **optimized for Molecular Devices Microplate Readers**. Click [here](#) for a list of recommended Microplate Readers.

If using a Molecular Devices' plate reader supported by SoftMax® Pro software, a preconfigured protocol for these CatchPoint SimpleStep ELISA Kits is available with all the protocol and analysis settings at [www.softmaxpro.org](http://www.softmaxpro.org)

The CatchPoint SimpleStep ELISA employs an affinity tag labeled capture antibody and a reporter conjugated detector antibody which immunocapture the sample analyte in solution. This entire complex (capture antibody/analyte/detector antibody) is in turn immobilized via immunoaffinity of an anti-tag antibody coating the well. To perform the assay, samples or standards are added to the wells, followed by the antibody mix. After incubation, the wells are washed to remove unbound material. CatchPoint HRP Development Solution containing the Stoplight Red Substrate is added. During incubation, the substrate is catalyzed by HRP generating a fluorescent product. Signal is generated proportionally to the amount of bound analyte and the intensity is measured in a fluorescence plate reader at 530/570/590 nm Excitation/Cutoff/Emission.

#### 特記事項

CXCL9 is a small cytokine belonging to the CXC chemokine subfamily that lacks an ELR motif in front of the first cysteine. CXCL9, also known as MIG (Monokine Induced by Gamma Interferon) is a T-cell chemoattractant, which is induced by Interferon Gamma. This subfamily also includes Interferon Gamma Induced Protein 10 (IP-10 or CXCL10) and Interferon Inducible T-cell Alpha Chemoattractant (I-TAC or CXCL11) whose genes are located near the gene for CXCL9 on human chromosome 4, CXCL9, IP-10, and I-TAC all elicit their chemotactic functions by interacting with the G protein coupled chemokine receptor CXCR3 (GPR9 or CD183).

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

#### 試験プラットフォーム

Pre-coated microplate (12 x 8 well strips)

#### 製品の特性

##### 保存方法

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

内容	1 x 96 tests
100X Stoplight Red Substrate	1 x 120µl
10X Human CXCL9 Capture Antibody	1 x 600µl
10X Human CXCL9 Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
500X Hydrogen Peroxide (H2O2, 3%)	1 x 50µl
50X Cell Extraction Enhancer Solution (ab193971)	1 x 1ml
5X Cell Extraction Buffer PTR (ab193970)	1 x 10ml

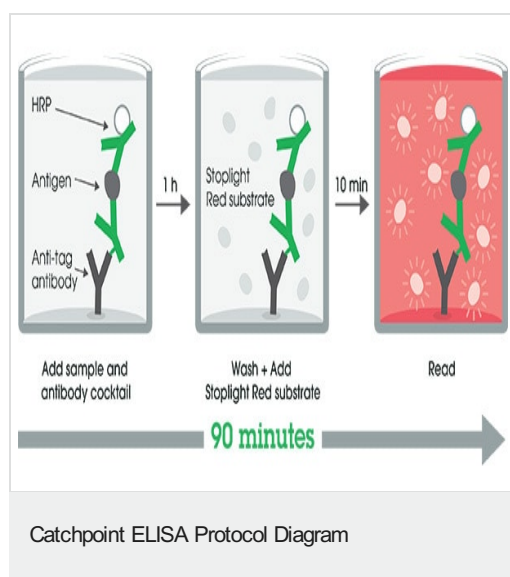
内容	1 x 96 tests
Antibody Diluent 4BI	1 x 6ml
Human CXCL9 Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated Black 96-Well Microplate	1 unit
Stoplight Red Substrate Buffer	1 x 12ml

**機能** Cytokine that affects the growth, movement, or activation state of cells that participate in immune and inflammatory response. Chemotactic for activated T-cells. Binds to CXCR3.

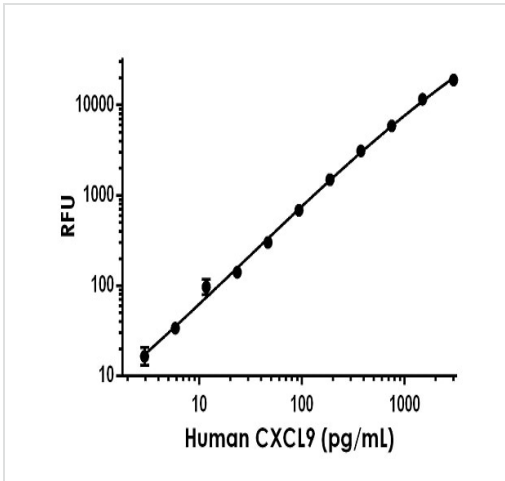
**配列類似性** Belongs to the intercrine alpha (chemokine CxC) family.

**細胞内局在** Secreted.

## 画像

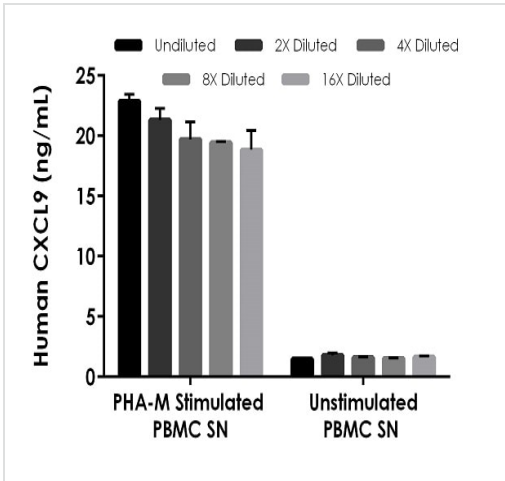


SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



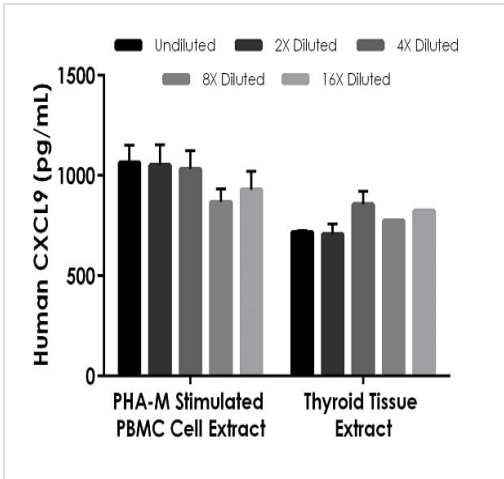
Example of human CXCL9 standard curve in Sample Diluent NS.

The CXCL9 standard curve was prepared as described in Section 10. Raw data generated on SpectraMax M4 Multi-Mode Microplate Reader is shown in the table. Background-subtracted data values (mean +/- SD) are graphed.



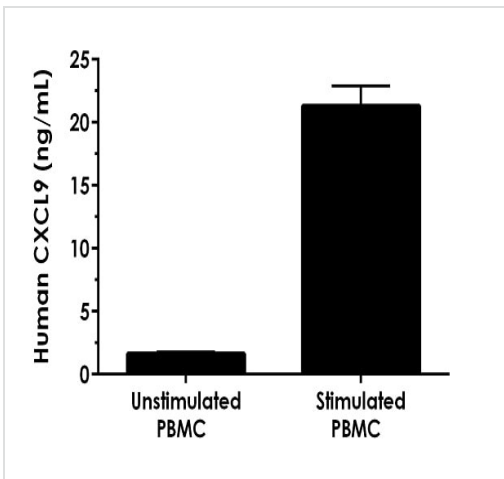
Interpolated concentrations of native CXCL9 in PHA-M stimulated and unstimulated human PBMC cell culture supernatant (2 days) samples.

The concentrations of CXCL9 were measured in duplicates, interpolated from the CXCL9 standard curves and corrected for sample dilution. Undiluted samples are as follows: PHA-M stimulated PBMC supernatant 2.5% and unstimulated PBMC supernatant 100% (neat). The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean CXCL9 concentration was determined to be 20.4 ng/mL in neat PHA-M stimulated PBMC supernatant and 1.62 ng/mL in neat unstimulated PBMC supernatant.



Interpolated concentrations of native CXCL9 in PHA-M stimulated human PBMC cell extract based on a 200 µg/mL extract load

The concentrations of CXCL9 were measured in duplicate and interpolated from the CXCL9 standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean CXCL9 concentration was determined to be 993.9 pg/mL in PHA-M stimulated PBMC cell extract and 824.9 pg/mL in thyroid tissue extract.



Comparison of CXCL9 in unstimulated and PHA-M stimulated human PBMC cell supernatants.

Human PBMC cells were cultured in the absence or presence of 1.5% PHA-M for 2 days. The concentrations of CXCL9 were measured in three different dilutions of the supernatant samples in duplicates and interpolated from the CXCL9 standard curve. The interpolated values are plotted (mean +/- SD, n=3). The mean CXCL9 concentration was determined to be 21.3 ng/mL in PHA-M stimulated PBMC cell supernatant, 1.6 ng/mL in unstimulated supernatants and undetectable in media (not shown).

Powered by  
recombinant antibodies



**Research with confidence**  
Consistent and reproducible results



**Long-term and scalable supply**  
Recombinant technology



**Success from the first experiment**  
Confirmed specificity



**Ethical standards compliant**  
Animal-free production

To learn more about the advantages of recombinant antibodies see [here](#).

Sandwich ELISA - Human CXCL9 ELISA  
Kit, Fluorescent (ab270891)

**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