

Product datasheet

Human IP-10 ELISA Kit, Fluorescent ab229410

CatchPoint® SimpleStep ELISA®

3 Images

Overview

Product name Human IP-10 ELISA Kit, Fluorescent

Detection method Fluorescent

Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Media	9			5.1%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Media	3			11.1%

Sample type Cell culture supernatant, Serum, Hep Plasma, EDTA Plasma, Cit plasma

Assay type Sandwich (quantitative)

Sensitivity 1.1 pg/ml

Range 1.6 pg/ml - 1600 pg/ml

Recovery	Sample specific recovery		
	Sample type	Average %	Range
	Serum	103	102% - 105%
	Cell culture media	96	95% - 98%
	Hep Plasma	86	82% - 88%
	EDTA Plasma	101	99% - 103%
	Cit plasma	96	93% - 98%

Assay time 1h 30m

Assay duration One step assay

**Species reactivity****Reacts with:** Human**Product overview**

IP-10 (CXCL10) *in vitro* CatchPoint SimpleStep ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IP-10 (CXCL10) protein in human serum and plasma 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.

**Notes**

C-X-C motif chemokine 10 (CXCL10 or IP-10) is a small 10.8kD protein that is secreted by several cell types in response to interferon-gamma (IFN $\gamma$ ). These cell types include monocytes, endothelial cells and fibroblasts. Upon secretion, CXCL10 is cleaved into an 8.7kD biologically active protein to function in chemotaxis for T-cells, NK cells, monocytes/macrophages and dendritic cells. In addition, CXCL10 has antitumor activity through the inhibition of bone marrow colony formation and angiogenesis. CXCL10 elicits its effects by binding to the cell surface chemokine receptor 3 (CXCR3).

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.

**Platform**

Pre-coated microplate (12 x 8 well strips)

**Properties****Storage instructions**

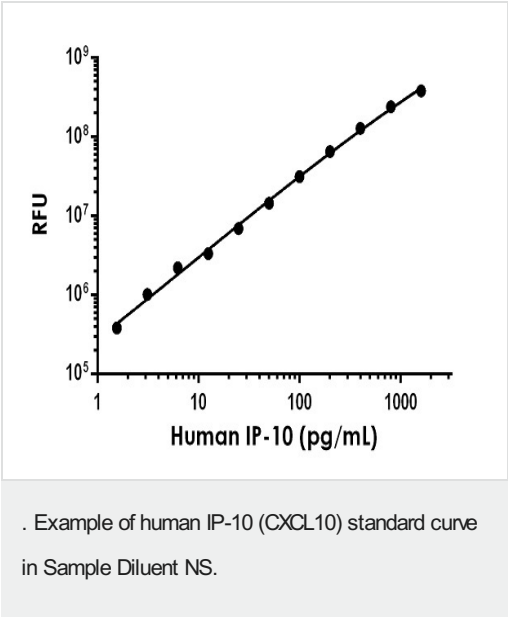
Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
100X Stoplight Red Substrate	1 x 120 $\mu$ l
10X Human IP-10 Capture Antibody	1 x 600 $\mu$ l
10X Human IP-10 Detector Antibody	1 x 600 $\mu$ l
10X Wash Buffer PT (ab206977)	1 x 20ml

Components	1 x 96 tests
500X Hydrogen Peroxide (H2O2, 3%)	1 x 50µl
Antibody Diluent CPI - HAMA Blocker (ab193969)	1 x 6ml
Human IP-10 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

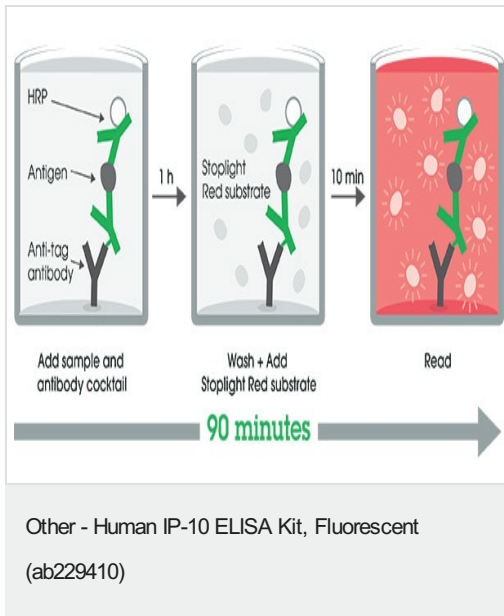
Function	Chemotactic for monocytes and T-lymphocytes. Binds to CXCR3.
Sequence similarities	Belongs to the intercrine alpha (chemokine CxC) family.
Post-translational modifications	CXCL10(1-73) is produced by proteolytic cleavage after secretion from keratinocytes.
Cellular localization	Secreted.

Images

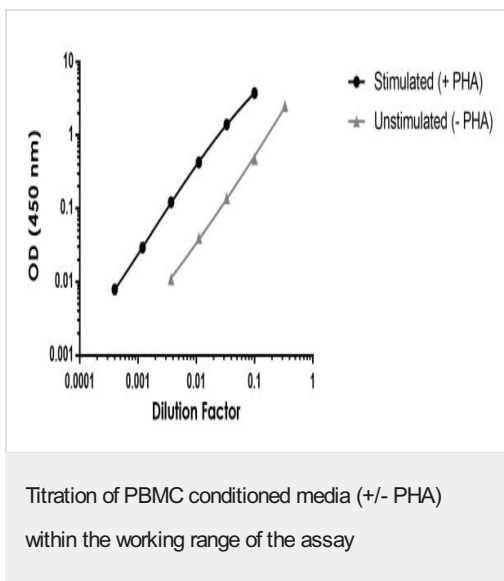


. Example of human IP-10 (CXCL10) standard curve in Sample Diluent NS.

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



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 from triplicate measurements are plotted.

**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.com/abpromise> or contact our technical team.

#### **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors