

Product datasheet

Human IP-10 ELISA Kit ab173194

SimpleStep ELISA[®]

★★★★★ 1 Abreviews 3 Images

Overview

Product name Human IP-10 ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
PBMC media	9			5.1%

Inter-assay

Sample	n	Mean	SD	CV%
PBMC media	3			11.1%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity 1.4 pg/ml

Range 1.4 pg/ml - 1000 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	103	102% - 105%
Cell culture media	96	95% - 98%
Heparin Plasma	86	82% - 88%
EDTA Plasma	101	99% - 103%
Citrate Plasma	96	93% - 98%

Assay time 1h 30m

Assay duration One step assay

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

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

The 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. TMB substrate is added and during incubation is catalyzed by HRP, generating blue coloration. This reaction is then stopped by addition of Stop Solution completing any color change from blue to yellow. Signal is generated proportionally to the amount of bound analyte and the intensity is measured at 450 nm. Optionally, instead of the endpoint reading, development of TMB can be recorded kinetically at 600 nm.

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 γ). 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).

Tested applications**Suitable for:** Sandwich ELISA**Platform**

Microplate

Properties**Storage instructions** Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
10X IP-10 Capture Antibody	1 x 600 μ l
10X IP-10 Detector Antibody	1 x 600 μ l
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent CPI - HAMA Blocker (ab193969)	1 x 6ml
IP-10 Human Lyophilized Recombinant Protein	2 x 1.25ng
Plate Seal	1 unit
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Substrate	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

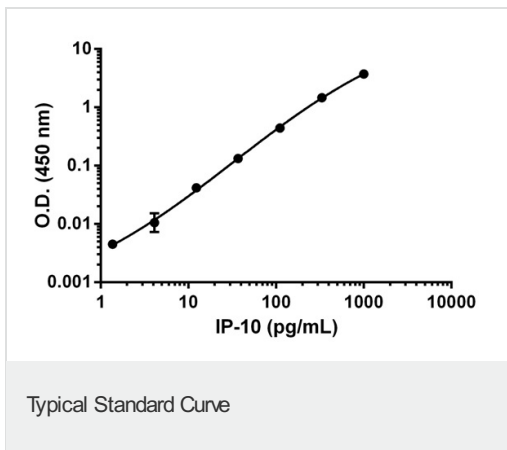
Secreted.

Applications

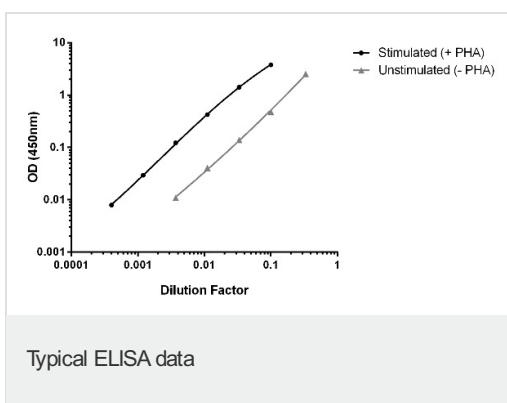
Our [Abpromise guarantee](#) covers the use of **ab173194** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

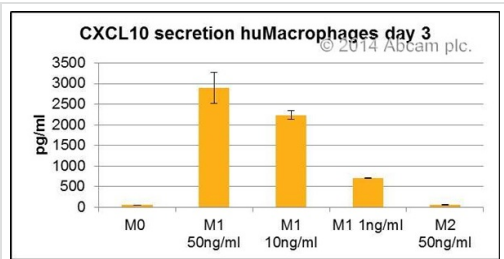
Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

Human IP-10 ELISA Kit images

Example IP-10/CXCL10 standard curve. Background-subtracted data values (mean +/- SD) are graphed. A new standard curve must be generated for each assay performed.



Titration of PBMC conditioned media (+/- PHA) within the working range of the assay. Background subtracted data from triplicate measurements are plotted.



Data shows specific secretion of IP-10 (CXCL10) by human macrophages differentiated in culture for 3 days in a dose response to M1 (MCSF + IFN γ).

Sandwich ELISA _ IP-10 (CXCL10) Human

SimpleStep ELISA™ Kit (ab173194)

This image is courtesy of an anonymous Abreview

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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