

## Product datasheet

# Mouse IP-10 ELISA Kit, Fluorescent ab272186

SimpleStep ELISA

[2 Images](#)

### Overview

**Product name** Mouse IP-10 ELISA Kit, Fluorescent

**Detection method** Fluorescent

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
Supernatant	8			8%

Inter-assay

Sample	n	Mean	SD	CV%
Supernatant	3			10.6%

**Sample type** Cell culture supernatant, Serum, Cit plasma

**Assay type** Sandwich (quantitative)

**Sensitivity** 50.32 pg/ml

**Range** 64.22 pg/ml - 32880 pg/ml

**Recovery**

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	104	93% - 115%
Serum	95	92% - 102%
Cit plasma	89	84% - 99%

**Assay time** 1h 30m

**Assay duration** One step assay

**Species reactivity** **Reacts with:** Mouse

**Product overview** IP-10 *in vitro* CatchPoint SimpleStep ELISA kit is designed for the quantitative measurement of IP-10 (CXCL10) protein in mouse serum, plasma citrate, and cell culture supernatant 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).

**Platform** Pre-coated microplate (12 x 8 well strips)

## Properties

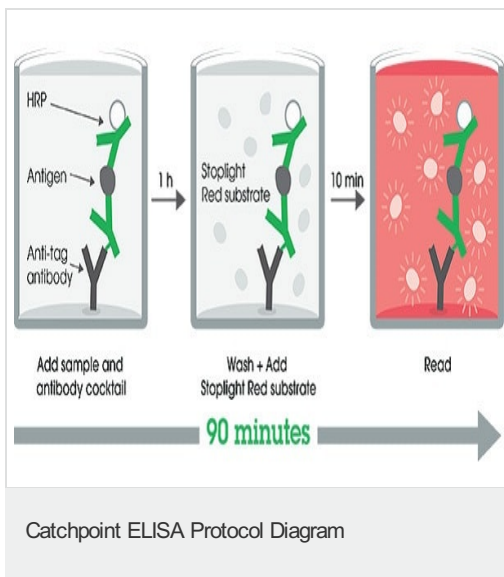
**Storage instructions** Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
100X Stoplight Red Substrate	1 x 120 $\mu$ l
10X Mouse IP-10 (CXCL10) Capture Antibody	1 x 600 $\mu$ l
10X Mouse IP-10 (CXCL10) Detector Antibody	1 x 600 $\mu$ l
10X Wash Buffer PT (ab206977)	1 x 20ml
500X Hydrogen Peroxide (H <sub>2</sub> O <sub>2</sub> , 3%)	1 x 50 $\mu$ l
Antibody Diluent 5BR	1 x 6ml
Mouse IP-10 (CXCL10) Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 12ml

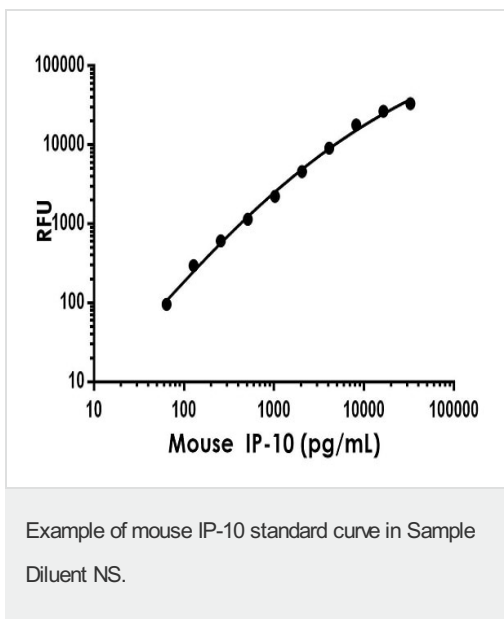
Components	1 x 96 tests
SimpleStep Pre-Coated Black 96-Well Microplate	1 unit
Stoplight Red Substrate Buffer	1 x 12ml

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

## Images



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.



The IP-10 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.

## **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