

Mouse IP-10 ELISA Kit, Fluorescent ab272186

SimpleStep ELISA

画像数 2

製品の概要

製品名 Mouse IP-10 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	8			8%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Supernatant	3			10.6%

サンプルの種類 Cell culture supernatant, Serum, Cit plasma

アッセイタイプ Sandwich (quantitative)

検出感度 50.32 pg/ml

検出範囲 64.22 pg/ml - 32880 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	104	93% - 115%
Serum	95	92% - 102%
Cit plasma	89	84% - 99%

全工程の試験時間 1h 30m

ステップ One step assay

種交差性 交差種: Mouse

製品の概要 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.

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

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).

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

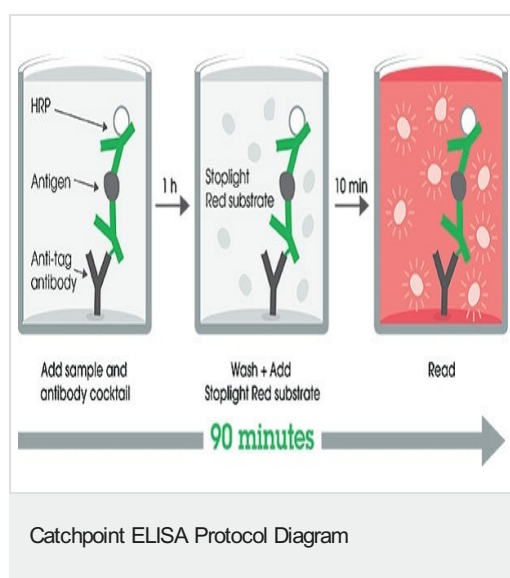
Store at +4°C. Please refer to protocols.

内容	1 x 96 tests
100X Stoplight Red Substrate	1 x 120 μ l
10X Mouse IP-10 (CXCL10) Capture Antibody	1 x 600 μ l
10X Mouse IP-10 (CXCL10) Detector Antibody	1 x 600 μ l
10X Wash Buffer PT (ab206977)	1 x 20ml
500X Hydrogen Peroxide (H ₂ O ₂ , 3%)	1 x 50 μ 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

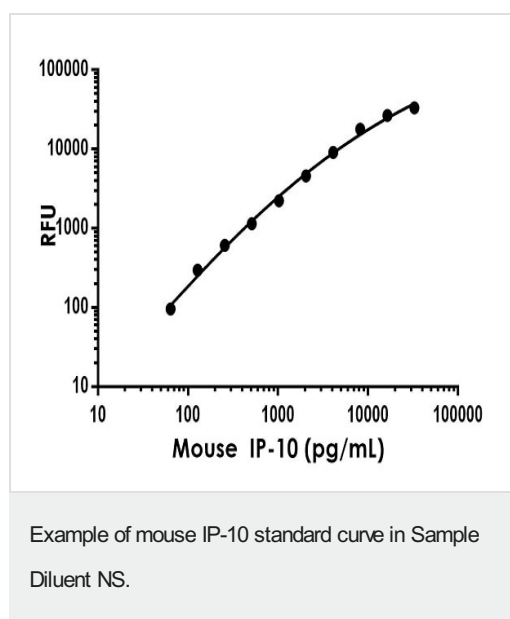
内容	1 x 96 tests
SimpleStep Pre-Coated Black 96-Well Microplate	1 unit
Stoplight Red Substrate Buffer	1 x 12ml

機能	Chemotactic for monocytes and T-lymphocytes. Binds to CXCR3.
配列類似性	Belongs to the intercrine alpha (chemokine CxC) family.
翻訳後修飾	CXCL10(1-73) is produced by proteolytic cleavage after secretion from keratinocytes.
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.



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.

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