

Human CSF-1-R ELISA Kit ab230940

リコンビナント SimpleStep ELISA

画像数 7

製品の概要

製品名 Human CSF-1-R ELISA Kit

検出方法 Colorimetric

再現性 Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
Human serum	9			6%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Human serum	9			5.6%

サンプルの種類 Serum, Cell culture extracts, Tissue Extracts, Hep Plasma, EDTA Plasma, Cit plasma

アッセイタイプ Sandwich (quantitative)

検出感度 20.1 pg/ml

検出範囲 78.1 pg/ml - 5000 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
Serum	105	104% - 107%
Cell culture extracts	102	101% - 103%
Tissue	108	106% - 110%
Tissue Culture Media	101	101% - 102%
Hep Plasma	108	105% - 111%
EDTA Plasma	108	106% - 111%

サンプルの種類	平均 %	測定範囲
Cit plasma	110	107% - 113%

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

非交差種: Cow, Macaque monkey

製品の概要

Human CSF-1-R ELISA Kit (ab230940) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of CSF-1-R protein in edta plasma, hep plasma, serum, tissue extracts, cit plasma, and cell culture extracts. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human CSF-1-R with 13.1 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate (**ab203359**) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

特記事項

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

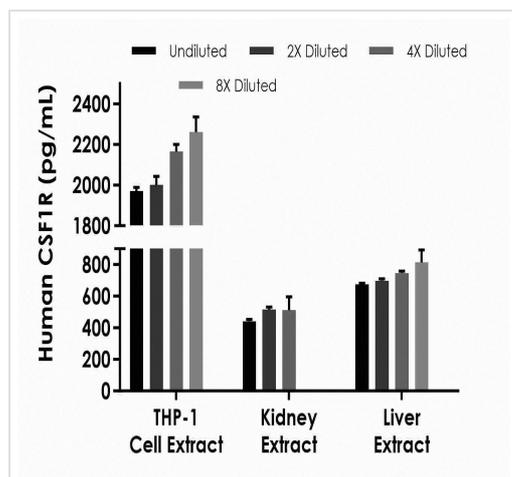
Store at +4°C. Please refer to protocols.

内容	1 x 96 tests
10X Human CSF-1-R Capture Antibody	1 x 600µl
10X Human CSF-1-R Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
50X Cell Extraction Enhancer Solution (ab193971)	1 x 1ml

内容	1 x 96 tests
5X Cell Extraction Buffer PTR (ab193970)	1 x 10ml
Antibody Diluent 4BR	1 x 6ml
Human CSF-1-R Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

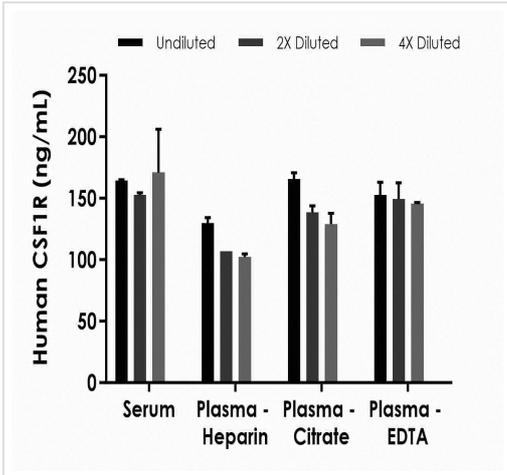
機能	Protein tyrosine-kinase transmembrane receptor for CSF1 and IL34.
組織特異性	Expressed in bone marrow and in differentiated blood mononuclear cells.
配列類似性	Belongs to the protein kinase superfamily. Tyr protein kinase family. CSF-1/PDGF receptor subfamily. Contains 5 Ig-like C2-type (immunoglobulin-like) domains. Contains 1 protein kinase domain.
細胞内局在	Membrane.

画像



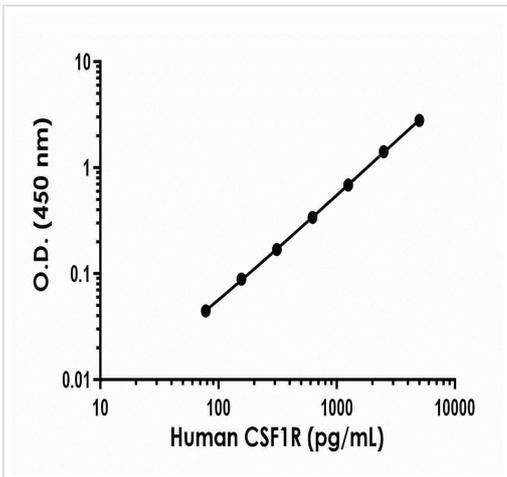
Interpolated concentrations of native CSF-1-R in human THP-1 cell extract, kidney extract, and liver extract

Interpolated concentrations of native CSF-1-R in human THP-1 cell extract, kidney extract, and liver extract. The concentrations of CSF-1-R were measured in duplicate and interpolated from the CSF-1-R standard curve and corrected for sample dilution. Undiluted samples are as follows: THP-1 cell extract 500 µg/mL, kidney extract 100 µg/mL, and liver extract 500 µg/mL. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean CSF-1-R concentration in the undiluted samples was determined to be 2,100 pg/mL in THP-1 cell extract, 489 pg/mL in kidney extract, and 733 pg/mL in liver extract.



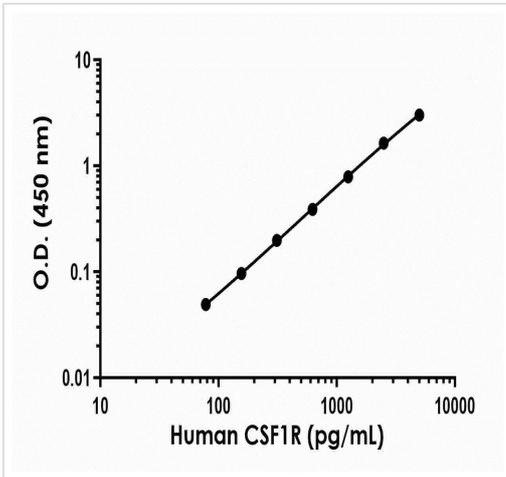
Interpolated concentrations of native CSF-1-R in human serum and plasma samples

Replace. Interpolated concentrations of native CSF-1-R in human serum and plasma samples. The concentrations of CSF-1-R were measured in duplicates, interpolated from the CSF-1-R standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 1:400, plasma (heparin) 1:200, plasma (citrate) 1:400 and plasma (EDTA) 1:400. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean CSF-1-R concentration was determined to be 163 ng/mL in neat serum, 113 ng/mL in neat plasma (heparin), 144 ng/mL in neat plasma (citrate), and 149 ng/mL in neat plasma (EDTA).



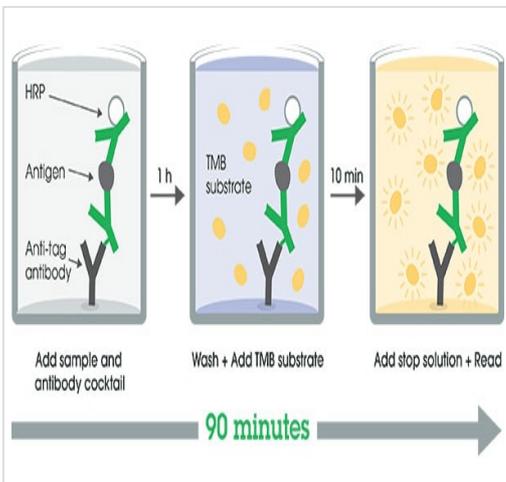
example of human CSF-1-R standard curve in Sample Diluent NS with Enhancer

The CSF-1-R standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.



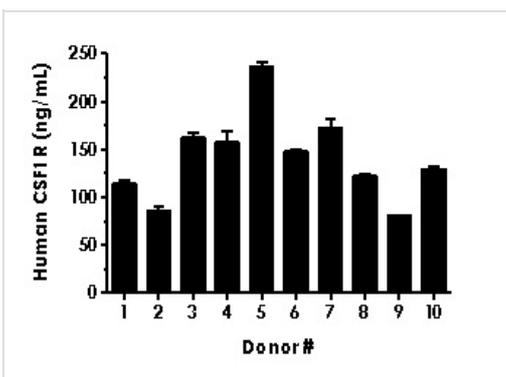
Example of human CSF-1-R standard curve in Sample Diluent NS

The CSF-1-R standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.



Other - Human CSF-1-R ELISA Kit (ab230940)

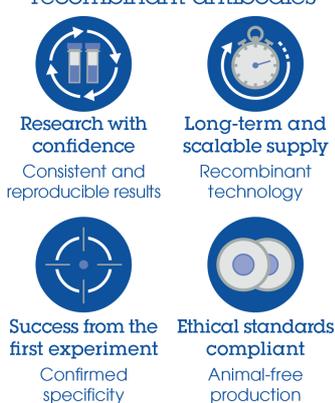
SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



Serum from ten individual healthy human female donors was measured in duplicate

Interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean CSF-1-R concentration was determined to be 141 ng/mL with a range of 81 – 240 ng/mL.

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

Sandwich ELISA - Human CSF-1-R ELISA Kit
(ab230940)

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

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