

Human Flt-3 ligand ELISA Kit ab254507

リコンビナント SimpleStep ELISA

画像数 8

製品の概要

製品名 Human Flt-3 ligand 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%
Serum	8			4.1%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Serum	3			4.2%

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

アッセイタイプ Sandwich (quantitative)

検出感度 2 pg/ml

検出範囲 4.7 pg/ml - 300 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	95	94% - 97%
Serum	99	93% - 103%
Hep Plasma	98	93% - 101%
EDTA Plasma	94	87% - 97%
Cit plasma	97	93% - 101%

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

製品の概要

Human Flt-3 ligand ELISA Kit (ab254507) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Flt-3 ligand protein in cell culture supernatant, cit plasma, edta plasma, hep plasma, and serum. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human Flt-3 ligand with 2 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

Flt-3 Ligand binds to and activates Flt-3. Flt-3 Ligand works synergistically with other CSFs and interleukins to stimulate growth and differentiation of early hematopoietic cells. The predominant form of Flt-3 Ligand is the biologically active transmembrane form which is proteolytically cleaved to generate a soluble form that is also biologically active. Alternative spliced secreted soluble variant of Flt-3 Ligand has also been identified.

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.

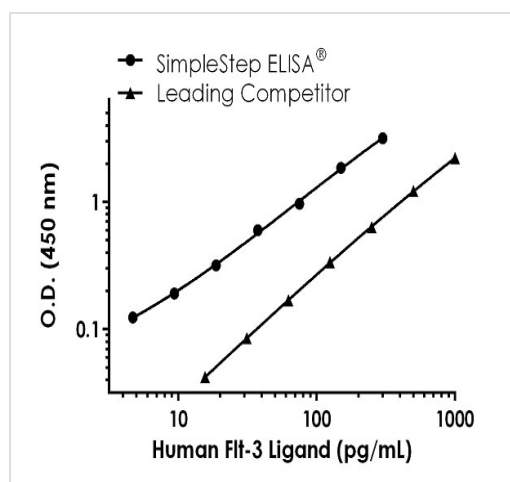
内容	1 x 96 tests	10 x 96 tests
10X Human Flt-3 ligand Capture Antibody	1 x 600µl	1 x 6000µl
10X Human Flt-3 ligand Detector Antibody	1 x 600µl	1 x 6000µl
10X Wash Buffer PT (ab206977)	1 x 20ml	1 x 200ml
Antibody Diluent 4BI	1 x 6ml	10 x 6ml
Human Flt-3 ligand Lyophilized Recombinant Protein	2 vials	2 x 10 vials
Plate Seals	1 unit	1 x 10 units
Sample Diluent NS (ab193972)	1 x 50ml	2 x 250ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit	1 x 10 units

内容	1 x 96 tests	10 x 96 tests
Stop Solution	1 x 12ml	1 x 120ml
TMB Development Solution	1 x 12ml	1 x 120ml

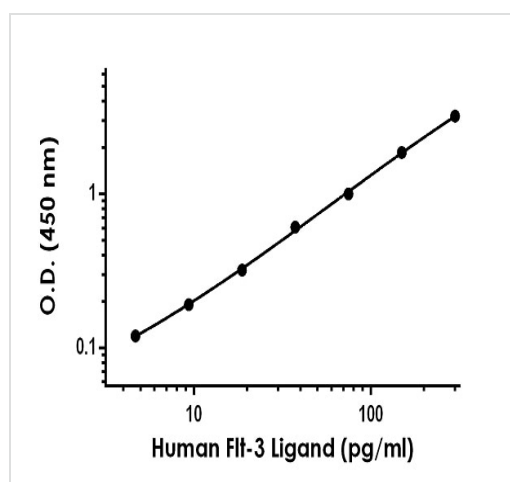
機能 Stimulates the proliferation of early hematopoietic cells by activating FLT3. Synergizes well with a number of other colony stimulating factors and interleukins.

細胞内局在 Secreted and Cell membrane.

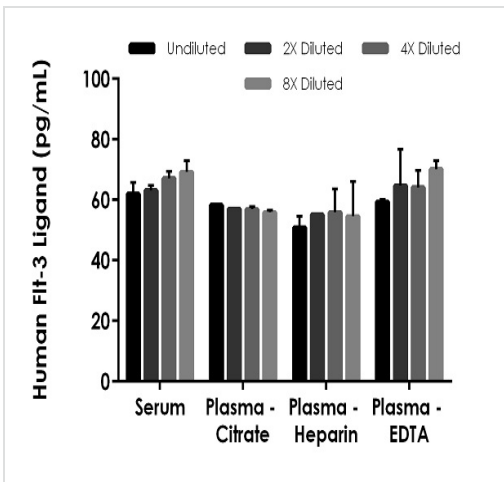
画像



Human Flt-3 Ligand standard curve comparison

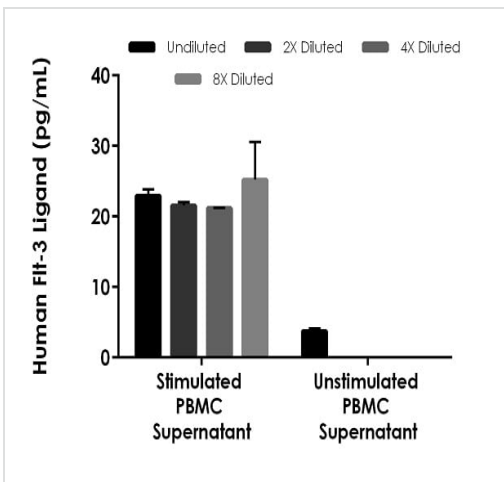


Example of human Flt-3 Ligand standard curve in Sample Diluent NS.



Interpolated concentrations of native Flt-3 Ligand in human serum and plasma samples.

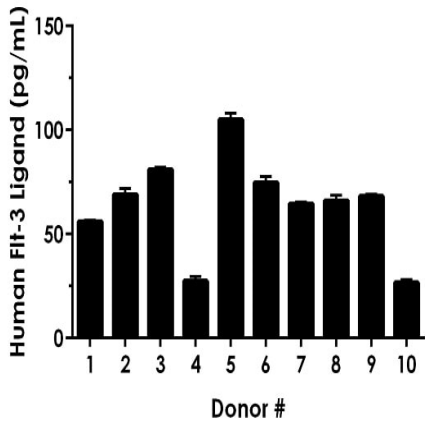
The concentrations of Flt-3 Ligand were measured in duplicates, interpolated from the Flt-3 Ligand standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 50%, plasma (citrate) 50%, plasma (heparin) 50%, and plasma (EDTA) 50%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Flt-3 Ligand concentration was determined to be 65.35 pg/mL in neat serum, 57.01 pg/mL in neat plasma (citrate), 54.11 pg/mL in neat plasma (heparin), and 64.57 pg/mL in neat plasma (EDTA).



Interpolated concentrations of native Flt-3 Ligand in human PBMC cell culture supernatant samples.

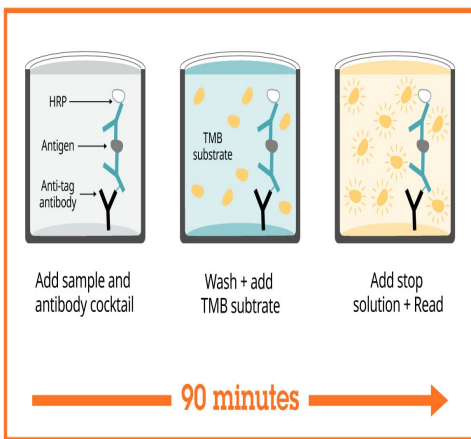
Human PBMC were cultured in the presence (stimulated) or absence (unstimulated) of 1.5% PHA-M.

Human peripheral blood mononuclear cells (PBMC) were cultured in the presence (stimulated) or absence (unstimulated) of 1.5% PHA-M for 36 hours. The concentrations of Flt-3 Ligand were measured in duplicates, interpolated from the Flt-3 Ligand standard curves and corrected for sample dilution. Undiluted samples are as follows: stimulated PBMC supernatant 100% and unstimulated PBMC supernatant 100%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Flt-3 Ligand concentration was determined to be 22.72 pg/mL in neat PHA-M stimulated PBMC cell culture supernatant, 3.742 pg/mL in neat unstimulated PBMC cell culture supernatant, and undetectable in naïve cell culture media (data not shown).



Interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Flt-3 Ligand concentration was determined to be 63.87 pg/mL with a range of 26.63 – 105.1 pg/mL.

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



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.

Sandwich ELISA - Human Flt-3 ligand ELISA Kit (ab254507)

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 Flt-3 ligand ELISA Kit
(ab254507)

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

Get more done with SimpleStep ELISA



Easy to use
Single-wash 90-minute protocol



Flexible
Matched antibody pairs available



Precision antibodies
High sensitivity, specificity and reproducibility



Scalable
Now in 10-pack and 384-well formats

Sandwich ELISA - Human Flt-3 ligand ELISA Kit
(ab254507)

To learn more about the advantages of SimpleStep ELISA[®] kits 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