

Mouse XCL1 ELISA Kit (Lymphotactin) ab269540

リコンビナント SimpleStep ELISA

画像数 7

製品の概要

製品名	Mouse XCL1 ELISA Kit (Lymphotactin)
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			4.5%

サンプル	N	平均値	SD	CV%
Serum	3			7.2%

サンプルの種類	Serum, Cell culture media, Hep Plasma, EDTA Plasma, Cit plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	6.26 pg/ml
検出範囲	18.75 pg/ml - 1200 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	116	110% - 123%
Cell culture media	99	94% - 105%
Hep Plasma	90	85% - 93%
EDTA Plasma	105	95% - 114%
Cit plasma	100	85% - 115%

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

Mouse XCL1 ELISA Kit (Lymphotactin) (ab269540) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of XCL1 (Lymphotactin) protein in cell culture media, edta plasma, hep plasma, serum, and cit plasma. It uses our proprietary SimpleStep ELISA® technology. Quantitate Mouse XCL1 (Lymphotactin) with 6.26 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

XCL1 (Lymphotactin) is a secreted glycosylated protein of C chemokine family. It plays a role in chemotactic activity of lymphocytes, but not monocytes or neutrophils. In the thymus, XCL1 mediates medullary accumulation of dendritic cells and contributes to T cell development which plays a role in establishment of self-tolerance. XCL2 is functionally related XCL1 homolog, these proteins share 98% amino acids sequence identity.

試験プラットフォーム

Pre-coated microplate (12 x 8 well strips)

製品の特性

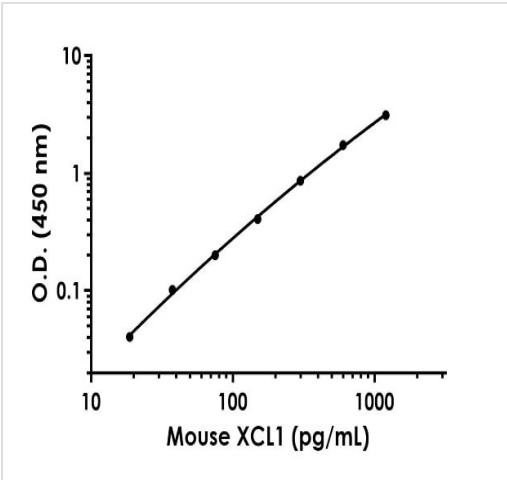
保存方法

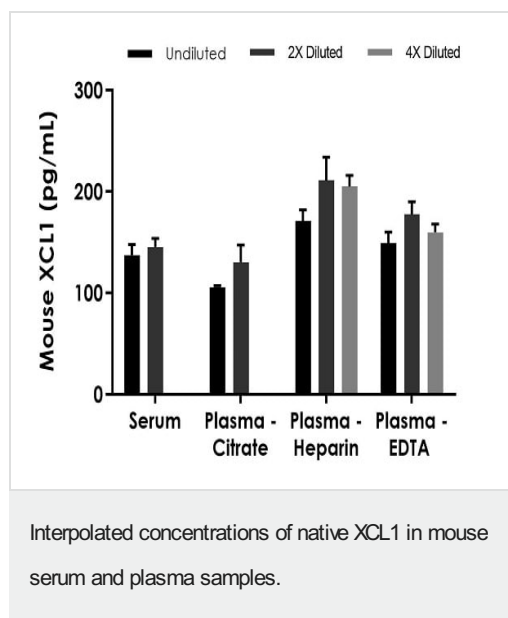
Store at +4°C. Please refer to protocols.

内容	1 x 96 tests
10X Mouse XCL1 (Lymphotactin) Capture Antibody	1 x 600µl
10X Mouse XCL1 (Lymphotactin) Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 4BR	1 x 6ml
Mouse XCL1 (Lymphotactin) Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 12ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit

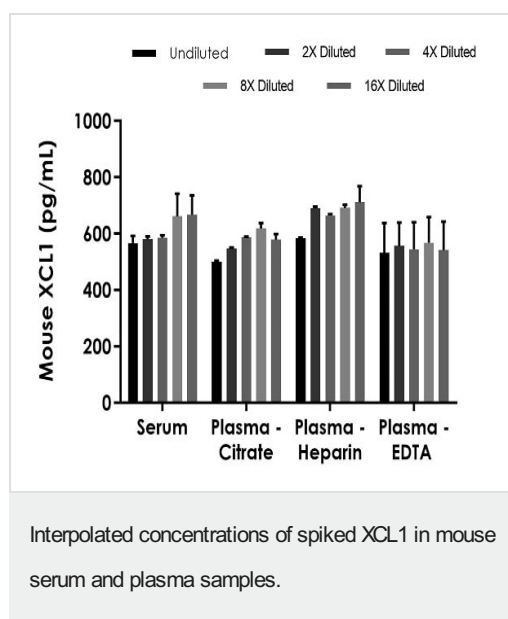
内容	1 x 96 tests
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

<p>関連性</p> <p>Human lymphotactin (Lptn)/XCL1, also named human SCM1 and ATAC, is a member of gamma or C subfamily of chemokines that lack two of the four invariant cysteine residues normally found in the CC and CXC chemokines and have an extended C-terminus. Human lymphotactin encodes a 114 amino acid precursor protein with a 21 amino acid predicted signal peptide. The gene for human and mouse lymphotactin has been mapped to chromosome 1. Human and mouse lymphotactin share approximately 60% amino acid sequence identity. The expression of lymphotactin is restricted to activated T cells such as activated CD8+ T cells and other class I MHC restricted T cells. Since lymphotactin is produced by lymphocytes and acts on lymphocytes, it is speculated that it is a messenger in T cell chemoattraction. Human lymphotactin is chemotactic for lymphocytes and NK cells. Its expression is abundant in some activated T cells such as activated CD8+ T cells and other class I MCH restricted T cells and is absent in CD4+ T cells. It has been shown that lymphotactin is a key regulator of lymphocyte motility and adhesion during acute allograft rejection. The specific receptor for lymphotactin is the orphan receptor XCR1 (GPR5). It has chemotactic activity for lymphocytes and natural killer cells but not for monocytes or neutrophils.</p>	
<p>細胞内局在</p>	<p>Secreted</p>

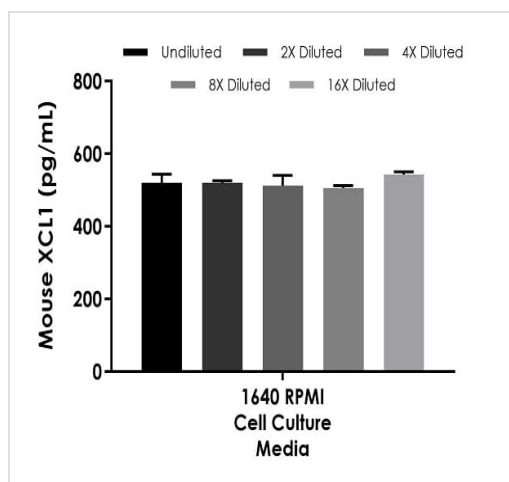
<p>画像</p>	
 <p>Example of mouse XCL1 standard curve in Sample Diluent NS.</p>	<p>The XCL1 standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.</p>



The concentrations of XCL1 were measured in duplicates, interpolated from the XCL1 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 \pm SD, n=2). The mean XCL1 concentration was determined to be 141.47 pg/mL in serum, 118.18 pg/mL in plasma (citrate), 195.80 pg/ml in plasma (heparin), and 162.36 pg/ml in plasma (EDTA).

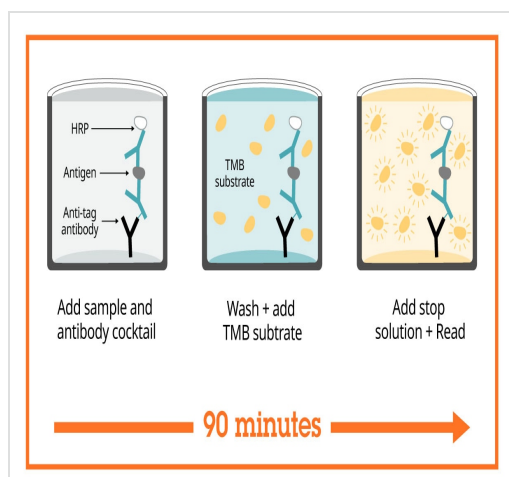


The concentrations of XCL1 were measured in duplicates, interpolated from the XCL1 standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 50%, plasma (citrate) 50%, plasma (heparin) 50%, and plasma (EDTA) 50%. The interpolated dilution factor corrected values are plotted (mean \pm SD, n=2).



Interpolated concentrations of spiked XCL1 in 1640 RPMI cell culture media samples containing 10% fetal bovine serum.

The concentrations of XCL1 were measured in duplicates, interpolated from the XCL1 standard curves and corrected for sample dilution. Undiluted samples are as follows: 1640 RPMI media 50%. The interpolated dilution factor corrected values are plotted (mean \pm SD, n=2).



Sandwich ELISA - Mouse XCL1 ELISA Kit
(Lymphotoxin) (ab269540)

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.

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 - Mouse XCL1 ELISA Kit
(Lymphotactin) (ab269540)

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 - Mouse XCL1 ELISA Kit
(Lymphotactin) (ab269540)

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