

Mouse CD30 ELISA Kit ab264614

リコンビナント SimpleStep ELISA

画像数 5

製品の概要

製品名 Mouse CD30 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%
Supernatant	8			3.5%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Supernatant	4			4.4%

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 79 pg/ml

検出範囲 133 pg/ml - 8500 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	106	86% - 119%
Serum	107	98% - 114%
Hep Plasma	96	93% - 108%
EDTA Plasma	99	90% - 106%
Cit plasma	95	87% - 102%

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 CD30 ELISA Kit (ab264614) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of CD30 protein in cell culture supernatant, cit plasma, edta plasma, hep plasma, and serum. It uses our proprietary SimpleStep ELISA® technology. Quantitate Mouse CD30 with 79 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

CD30, also known as Tumor necrosis factor receptor superfamily member 8, is a cell membrane receptor encoded by the gene *Tnfrsf8*. May play a role in the regulation of cellular growth and transformation of activated lymphoblasts.

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**

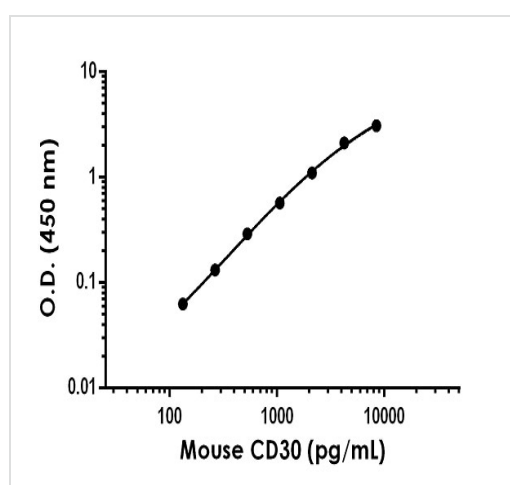
Store at +4°C. Please refer to protocols.

内容	1 x 96 tests
10X Mouse CD30 Capture Antibody	1 x 600µl
10x Mouse CD30 Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 4BR	1 x 6ml
Mouse CD30 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
Stop Solution	1 x 12ml

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

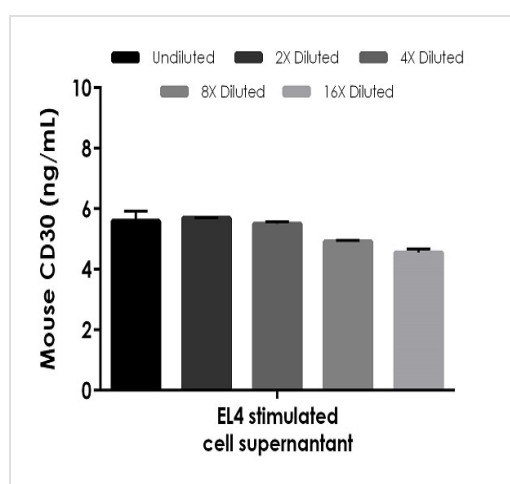
機能	Receptor for TNFSF8/CD30L. May play a role in the regulation of cellular growth and transformation of activated lymphoblasts. Regulates gene expression through activation of NF-kappa-B.
配列類似性	Contains 6 TNFR-Cys repeats.
翻訳後修飾	Phosphorylated on serine and tyrosine residues.
細胞内局在	Cytoplasm and Cell membrane.

画像



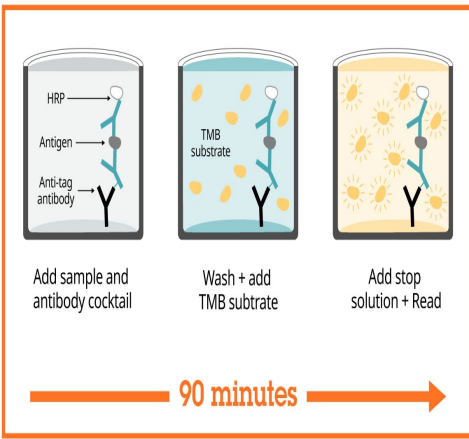
Example of mouse CD30 standard curve in Sample Diluent NS.

The CD30 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.



Interpolated concentrations of native CD30 in mouse cell culture supernatant samples.

The concentrations of CD30 were measured in duplicates, interpolated from the CD30 standard curves and corrected for sample dilution. Undiluted samples are as follows: Stimulated EL4 supernatant 100%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean CD30 concentration was determined to be 5.3 ng/mL in stimulated EL4 supernatant.



Sandwich ELISA - Mouse CD30 ELISA Kit
(ab264614)

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 CD30 ELISA Kit
(ab264614)

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 CD30 ELISA Kit
(ab264614)

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