

Mouse CD40L ELISA Kit ab119517

4 References [画像数 1](#)

製品の概要

製品名 Mouse CD40L 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%
Overall	8			6.5%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Overall	8			11.1%

サンプルの種類 Cell culture supernatant, Serum

アッセイタイプ Sandwich (quantitative)

検出感度 0.14 ng/ml

検出範囲 0.31 ng/ml - 20 ng/ml

添加回収試験 94 %

ステップ Multiple steps standard assay

種交差性 交差種: Mouse

製品の概要 Abcam's CD40L Mouse in vitro ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for accurate quantitative measurement of Mouse CD40L concentrations in cell culture supernatant and serum.

CD40L specific antibodies have been precoated onto 96-well plates. Standards and test samples are added to the wells along with a biotin-conjugated CD40L detection antibody then incubated at room temperature. Following washing, a Streptavidin-HRP conjugate is added to each well, incubated at room temperature and washed. TMB is added and then catalyzed by HRP to produce a blue product that changes to yellow after the addition of acidic stop solution. The density of yellow coloration is directly proportional to the amount of CD40L captured on the plate.

試験プラットフォーム Microplate

保存方法 Store at +4°C. Please refer to protocols.

内容	1 x 96 tests
20X Assay Buffer Concentrate	1 x 5ml
20X Wash Buffer Concentrate	1 x 50ml
Adhesive Films	4 units
Biotin-Conjugate anti-mouse CD40L monoclonal antibody	1 x 70µl
Microplate coated with monoclonal antibody to mouse CD40L (12 x 8 wells)	1 unit
Mouse CD40L Standard lyophilized (40 ng /mL upon reconstitution)	2 vials
Sample Diluent	1 x 12ml
Stop Solution (1M Phosphoric acid)	1 x 15ml
Streptavidin-HRP	1 x 150µl
TMB Substrate Solution	1 x 15ml

機能 Mediates B-cell proliferation in the absence of co-stimulus as well as IgE production in the presence of IL-4. Involved in immunoglobulin class switching.
Release of soluble CD40L from platelets is partially regulated by GP IIb/IIIa, actin polymerization, and an matrix metalloproteinases (MMP) inhibitor-sensitive pathway.

組織特異性 Specifically expressed on activated CD4+ T-lymphocytes.

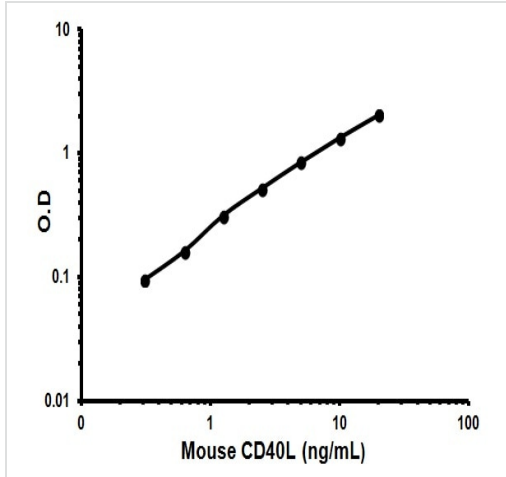
関連疾患 Defects in CD40LG are the cause of X-linked immunodeficiency with hyper-IgM type 1 (HIGM1) [MIM:308230]; also known as X-linked hyper IgM syndrome (XHIM). HIGM1 is an immunoglobulin isotype switch defect characterized by elevated concentrations of serum IgM and decreased amounts of all other isotypes. Affected males present at an early age (usually within the first year of life) recurrent bacterial and opportunistic infections, including Pneumocystis carinii pneumonia and intractable diarrhea due to cryptosporidium infection. Despite substitution treatment with intravenous immunoglobulin, the overall prognosis is rather poor, with a death rate of about 10% before adolescence.

配列類似性 Belongs to the tumor necrosis factor family.

翻訳後修飾 The soluble form derives from the membrane form by proteolytic processing.
N-linked glycan is a mixture of high mannose and complex type. Glycan structure does not influence binding affinity to CD40.
Not O-glycosylated.

細胞内局在 Secreted and Cell membrane.

画像



Representative Standard Curve using ab119517.

Typical Standard Curve

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