

Human CD40L ELISA Kit ab99991

3 References [画像数 2](#)

製品の概要

製品名	Human CD40L ELISA Kit
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 6 pg/ml
検出範囲	8.23 pg/ml - 6000 pg/ml
添加回収試験	95 %

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	96.35	86% - 105%
Serum	93.57	83% - 103%
Plasma	94.63	84% - 105%

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要

Human CD40L ELISA Kit (ab99991) is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human CD40L in serum, plasma (EDTA or heparin), and cell culture supernatants.

This assay employs an antibody specific for Human CD40L coated on a 96-well plate. Standards and samples are pipetted into the wells and CD40L present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human CD40L antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of CD40L bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Get results in 90 minutes with Human CD40 Ligand ELISA Kit (TNFSF5) ([ab196268](#)) from our SimpleStep ELISA[®] range.

特記事項 Optimization may be required with urine samples.

試験プラットフォーム Microplate

製品の特性

保存方法 Store at -20°C. Please refer to protocols.

内容	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
600X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated anti-Human CD40L	2 vials
CD40L Microplate (12 x 8 well strips)	1 unit
Recombinant Human CD40L Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

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.

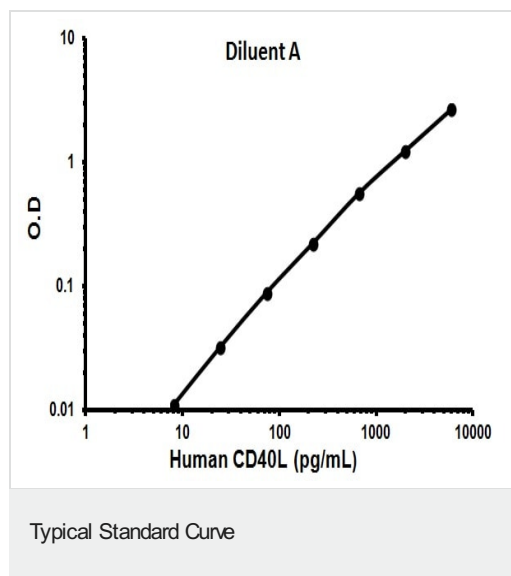
組織特異性 Specifically expressed on activated CD4+ T-lymphocytes.

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.

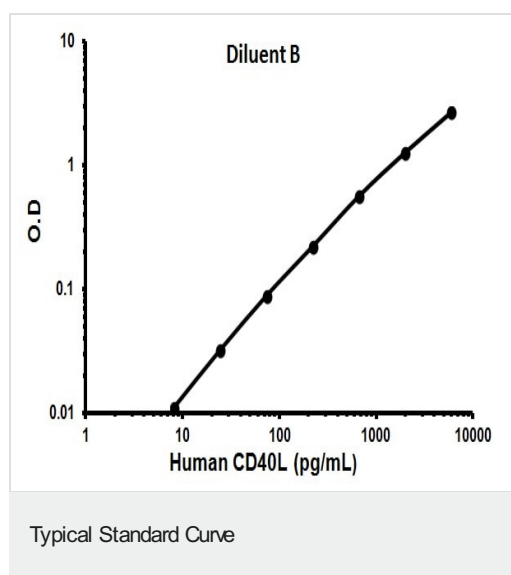
配列類似性 Belongs to the tumor necrosis factor family.

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.



Representative Standard Curve using ab99991.



Representative Standard Curve using ab99991.

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