

Human CD40L ELISA Kit ab99991

3 References 2 图像

概述

产品名称	人CD40L ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 6 pg/ml
范围	8.23 pg/ml - 6000 pg/ml
回收率	95 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	96.35	86% - 105%
Serum	93.57	83% - 103%
Plasma	94.63	84% - 105%

实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	<p>Human CD40L ELISA Kit (ab99991) is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human CD40L in serum, plasma (EDTA or heparin), and cell culture supernatants.</p> <p>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.</p> <p>Get results in 90 minutes with Human CD40 Ligand ELISA Kit (TNFSF5) (<a href="#">ab196268</a>) from our SimpleStep ELISA® range.</p>

说明 Optimization may be required with urine samples.

平台 Microplate

性能

存放说明 Store at -20°C. Please refer to protocols.

组 件	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

功能 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.

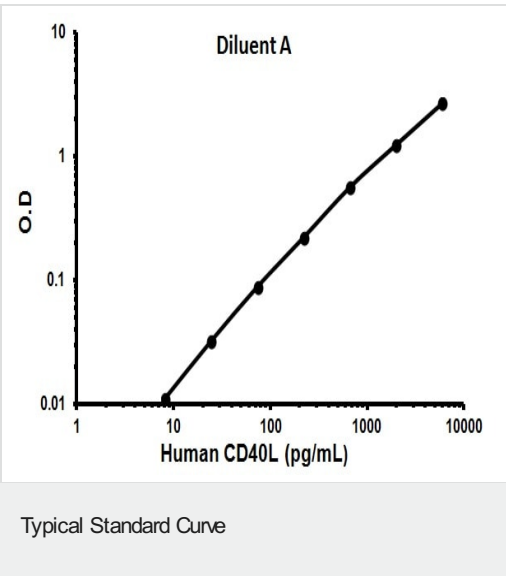
组织特异性 Specifically expressed on activated CD4+ T-lymphocytes.

疾病相关 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.

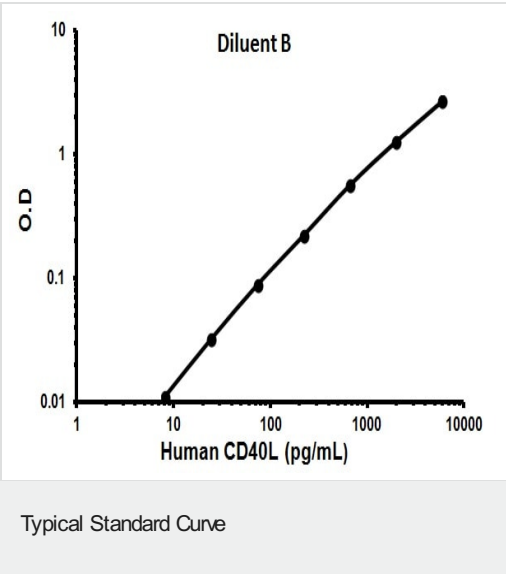
序列相似性 Belongs to the tumor necrosis factor family.

翻译后修饰 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.

细胞定位 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.cn/abpromise> or contact our technical team.

## Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors