

## Product datasheet

# Human CD40 ELISA Kit ab119516

### 画像数 1

#### 製品の概要

**製品名** Human CD40 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			5.5%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Overall	8			7%

**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** 1.3 pg/ml

**検出範囲** 7.8 pg/ml - 500 pg/ml

**添加回収試験** 94 %

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	94	81% - 110%

**ステップ** Multiple steps standard assay

**種交差性** 交差種: Human

**製品の概要** Abcam's CD40 Human in vitro ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for accurate quantitative measurement of Human CD40 concentrations in cell culture supernatant, serum, plasma (EDTA, heparin) and amniotic fluid.

CD40 specific antibodies have been precoated onto 96-well plates. Standards and test samples are added to the wells along with a biotinylated CD40 detection antibody then incubated at room temperature and washed. A Streptavidin-HRP conjugate is then added to each well, incubated at room temperature and washed. TMB is added and catalyzed by HRP to

produce a blue color product that changes into yellow after addition of acidic stop solution. The density of yellow coloration is directly proportional to the amount of CD40 sample captured in plate.

#### アプリケーション

**適用あり:** Sandwich ELISA

#### 試験プラットフォーム

Microplate

#### 製品の特性

#### 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-Human CD40 monoclonal antibody	1 x 70µl
CD40 Standard Lyophilized (1000 pg/mL on reconstitution)	2 vials
Microplate coated with anti-Human CD40 monoclonal antibody (12 x 8 wells)	1 unit
Sample Diluent	1 x 12ml
Stop Solution (1M Phosphoric acid)	1 x 15ml
Streptavidin-HRP	1 x 150µl
TMB Substrate Solution	1 x 15ml

#### 関連性

CD40 (48 to 50 kDa) is a transmembrane glycoprotein mainly expressed on the surface of B cells and also expressed on monocytes, dendritic cells, and thymic epithelium. CD40 is a member of the tumor necrosis factor (TNF) receptor superfamily, which includes the low affinity nerve growth factor (NGF) receptor and CD95/Fas. CD40 is the receptor for CD40 ligand. CD40 ligand (CD40L, CD154, gp39, and TRAM) belongs to the TNF gene family and is expressed more widely than CD40 predominantly on activated CD4+ T cells. Following interaction with CD40 ligand, CD40 mediates a number of major immunoregulatory functions, central to the control of thymus dependent humoral immunity and may be critical in the development of cell mediated immune responses. Other biological actions include B cell homotypic adhesion, proliferation, immunoglobulin isotype switch, and secretion. Activation of CD40 has also been shown to inhibit the growth of certain B cell lymphomas and to induce the death of transformed cells of mesenchymal or epithelial origin.

#### アプリケーション

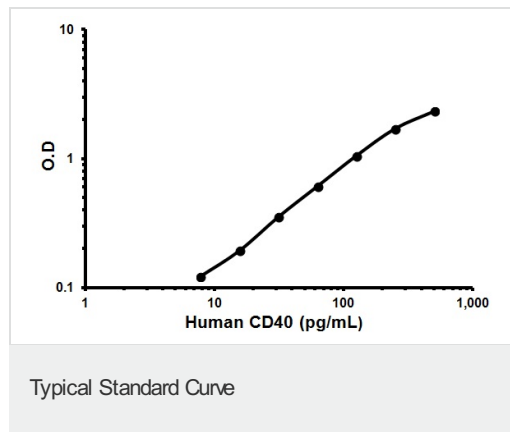
Our [Abpromise guarantee](#) covers the use of **ab119516** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Sandwich ELISA

Use at an assay dependent concentration.

## 画像



Representative Standard Curve using  
ab119516.

**Please note:** All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

### 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 <http://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