

Human VCAM1 ELISA Kit ab100661

3 References 画像数 2

製品の概要

|         |   |
|---------|---|
| 製品名     | Human VCAM1 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    | < 0.3 ng/ml                             |
| 検出範囲    | 0.247 ng/ml - 60 ng/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 | 98.76 | 85% - 104% |
| Serum                    | 94.42 | 83% - 103% |
| Plasma                   | 93.58 | 82% - 102% |

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 Abcam's VCAM1 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human VCAM1 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Human VCAM1 coated on a 96-well plate. Standards and samples are pipetted into the wells and VCAM1 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human VCAM1 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 VCAM1 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

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     |
| 320X HRP-Streptavidin Concentrate              | 1 x 200µl    |
| 5X Assay Diluent B                             | 1 x 15ml     |
| Assay Diluent A                                | 1 x 30ml     |
| Biotinylated anti-Human VCAM1                  | 2 vials      |
| Recombinant Human VCAM1 Standard (lyophilized) | 2 vials      |
| Stop Solution                                  | 1 x 8ml      |
| TMB One-Step Substrate Reagent                 | 1 x 12ml     |
| VCAM1 Microplate (12 x 8 wells)                | 1 unit       |

### 機能

Important in cell-cell recognition. Appears to function in leukocyte-endothelial cell adhesion. Interacts with the beta-1 integrin VLA4 on leukocytes, and mediates both adhesion and signal transduction. The VCAM1/VLA4 interaction may play a pathophysiologic role both in immune responses and in leukocyte emigration to sites of inflammation.

### 組織特異性

Expressed on inflamed vascular endothelium, as well as on macrophage-like and dendritic cell types in both normal and inflamed tissue.

### 配列類似性

Contains 7 Ig-like C2-type (immunoglobulin-like) domains.

### ドメイン

Either the first or the fourth Ig-like C2-type domain is required for VLA4-dependent cell adhesion.

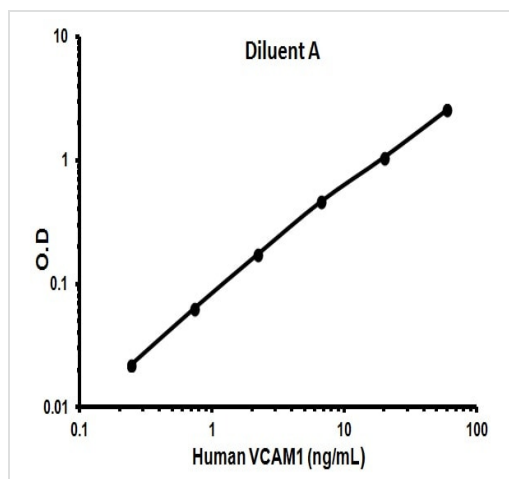
### 翻訳後修飾

Sialoglycoprotein.

### 細胞内局在

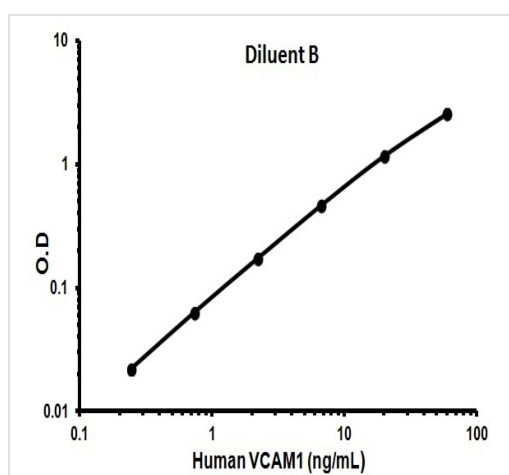
Membrane.

## 画像



Representative standard curve using ab100661

Typical standard curve



Representative standard curve using ab100661

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