

Mouse VCAM1 ELISA Kit ab100750

5 References 2 Images

Overview

Product name	Mouse VCAM1 ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 1.5 pg/ml
Range	2.05 pg/ml - 500 pg/ml
Recovery	96 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	97.36	84% - 107%
Serum	96.43	82% - 105%
Plasma	94.64	81% - 104%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview Abcam's VCAM1 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of mouse VCAM1 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Mouse 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-Mouse 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.

Platform Microplate

Properties

**Storage instructions**

Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
120X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent A	2 x 30ml
Biotinylated anti-Mouse VCAM1	2 vials
Recombinant Mouse VCAM1 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml
VCAM1 Microplate (12 strips x 8 wells)	1 unit

**Function**

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.

**Tissue specificity**

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

**Sequence similarities**

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

**Domain**

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

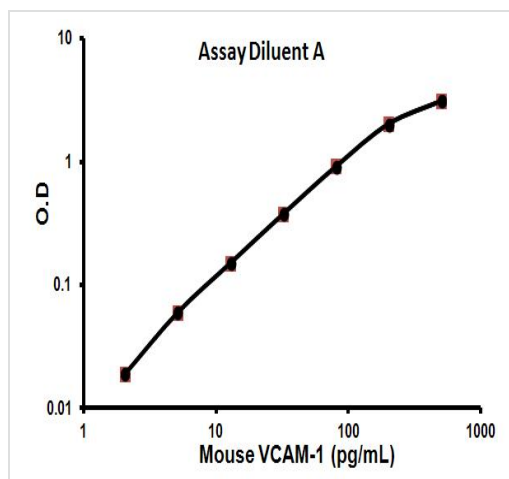
**Post-translational modifications**

Sialoglycoprotein.

**Cellular localization**

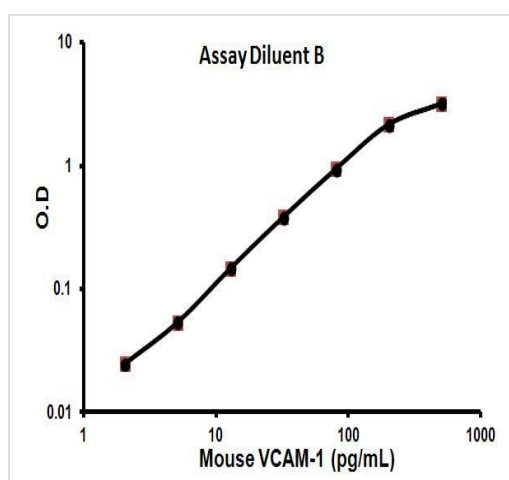
Membrane.

**Images**



Representative Standard Curve using ab100750

Representative Standard Curve



Representative Standard Curve using ab100750

Representative 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.com/abpromise> or contact our technical team.

### Terms and conditions

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