

### Human VCAM-1 ELISA Set (without plates) ab47355

[3 References](#) [1 Image](#)

#### Overview

<b>Product name</b>	Human VCAM-1 ELISA Set (without plates)
<b>Sample type</b>	Cell culture supernatant, Serum, Plasma
<b>Assay type</b>	Sandwich (quantitative)
<b>Sensitivity</b>	< 0.6 ng/ml
<b>Range</b>	1.56 ng/ml - 50 ng/ml
<b>Assay time</b>	1h 45m
<b>Assay duration</b>	Multiple steps standard assay
<b>Species reactivity</b>	<b>Reacts with:</b> Human
<b>Product overview</b>	Abcam's VCAM-1 Human <i>in vitro</i> ELISA Set is designed for the quantitative measurement of VCAM-1 in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.

This ELISA Set is intended to measure human VCAM-1 in cell -culture supernatants, as well as in serum and plasma samples. The enclosed procedures provide general guidelines for ELISA preparation. This procedure may need some optimization depending on the plastic, matrix, standard range.

The VCAM-1 ELISA Set includes both the capture and the detection antibodies as well as Horse-Radish peroxidase conjugated Streptavidin (HRP-Strep) and TMB which are the reagents used in the colour development step.

Recognizes both natural and recombinant human VCAM1

<b>Tested applications</b>	<b>Suitable for:</b> Sandwich ELISA
<b>Platform</b>	Reagents

#### Properties

**Storage instructions** Store at +4°C. Please refer to protocols.

Components	10 x 96 tests	15 x 96 tests	5 x 96 tests
Chromogen TMB Substrate Solution	4 vials	6 vials	2 vials

Components	10 x 96 tests	15 x 96 tests	5 x 96 tests
Detection biotinylated anti-VCAM-1 antibody	2 x 0.55ml	3 x 0.55ml	1 x 0.55ml
Streptavidin-HRP	2 vials	3 vials	1 vial
VCAM-1 Capture Antibody	2 x 0.5ml	3 x 0.5ml	1 x 0.5ml
VCAM1 standard	10 vials	15 vials	5 vials

**Relevance**

VCAM1 is important in cell-cell recognition. Appears to function in leukocyte-endothelial cell adhesion. Interacts with the integrins alpha4 beta1 (beta 1 integrin VLA4) and alpha4 beta7 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. VCAM1 is also expressed by several non endothelial cell types including some macrophages, follicular dendritic cells and bone marrow, stromal cells.

**Cellular localization**

Cell Membrane; single-pass type I membrane protein

**Applications**

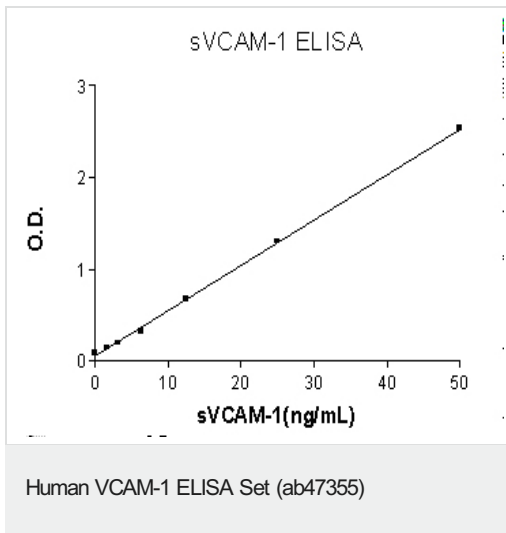
**The Abpromise guarantee**

Our **Abpromise guarantee** covers the use of ab47355 in the following tested applications.

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

Application	Abreviews	Notes
<b>Sandwich ELISA</b>		Use at an assay dependent concentration.

**Images**



VCAM-1 Standard curve ranging from 1.56 to 50 ng/mL

**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