

Human VCAM1 ELISA Kit (CD106) ab46118

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

Overview

Product name Human VCAM1 ELISA Kit (CD106)

Detection method Colorimetric

Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	A	8	48.79ng/ml	0.221	0.45%
	B	8	5.81ng/ml	0.132	2.27%
	Inter-assay				
	Sample	n	Mean	SD	CV%

	Sample	n	Mean	SD	CV%
	A	28	49.69ng/ml	0.72	1.44%
	B	28	5.95ng/ml	0.354	5.94%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity < 0.6 ng/ml

Range 1.56 ng/ml - 50 ng/ml

Assay time 1h 45m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's Human VCAM1 (CD106) *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of VCAM1 in cell culture supernatants, buffered solutions, serum, plasma, and other body fluids.

A monoclonal antibody specific for VCAM1 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known VCAM1 concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for VCAM1 are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated

and washed. A TMB substrate solution is added which acts on the bound enzyme to induce a colored reaction product. The intensity of this colored product is directly proportional to the concentration of VCAM1 present in the samples.

This kit will recognize both endogenous and recombinant Human VCAM1.

Get results in 90 minutes with Human VCAM1 ELISA Kit (CD106) (**ab187393**) from our SimpleStep ELISA® range.

Platform Microplate

Properties

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

Components	Identifier	1 x 96 tests	2 x 96 tests	1 x 96 tests
10X Standard Diluent Buffer	Black	1 x 15ml	1 x 25ml	1 x 15ml
200X Wash Buffer	White	1 x 10ml	2 x 10ml	1 x 10ml
Biotinylated Antibody Diluent	Red	1 x 7.5ml	1 x 13ml	1 x 7.5ml
Biotinylated anti-VCAM1	Red	1 x 400µl	2 x 400µl	1 x 400µl
Chromogen TMB Substrate Solution		1 x 11ml	1 x 24ml	1 x 11ml
Control	Silver	2 vials	4 vials	2 vials
HRP Diluent	Red	1 x 12ml	1 x 23ml	1 x 12ml
Stop Reagent	Black	1 x 11ml	2 x 11ml	1 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl	2 x 5µl
VCAM1 Microplate (12 x 8 well strips)		1 unit	2 units	1 unit
VCAM1 Standard (lyophilized)	Yellow	2 vials	4 vials	2 vials

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.

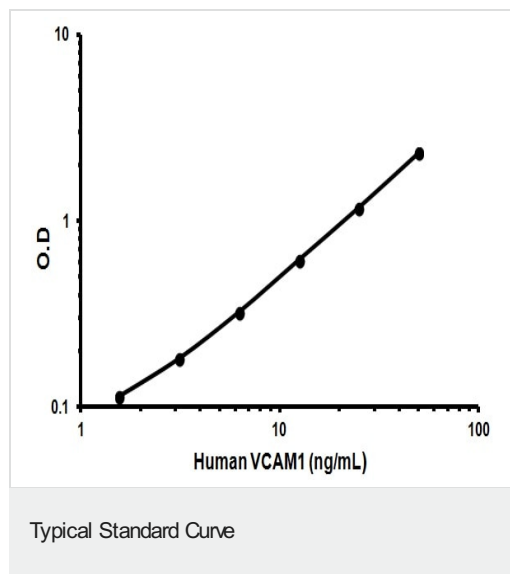
Post-translational Sialoglycoprotein.

modifications

Cellular localization

Membrane.

Images



Representative Standard Curve using ab46118.

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