

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

3 References 1 图像

概述

产品名称	人VCAM-1 ELISA试剂盒(without plates)
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 0.6 ng/ml
范围	1.56 ng/ml - 50 ng/ml
检测时间	1h 45m
实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	<p>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.</p> <p>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.</p> <p>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.</p> <p>Recognizes both natural and recombinant human VCAM1</p>
经测试应用	适用于: Sandwich ELISA
平台	Reagents

性能

存放说明 Store at +4°C. Please refer to protocols.

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

组件	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

相关性	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.
细胞定位	Cell Membrane; single-pass type I membrane protein

应用

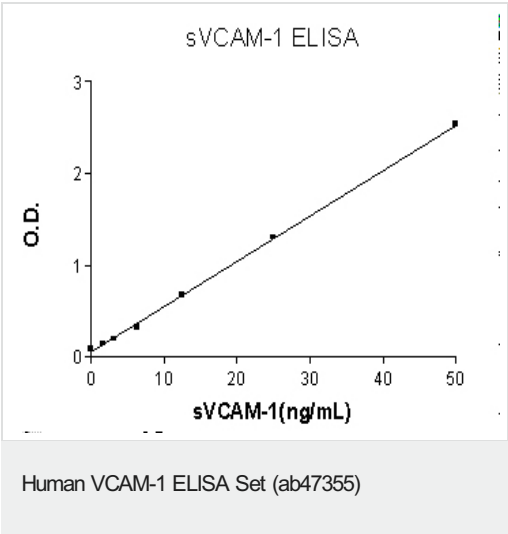
The Abpromise guarantee

Abpromise™承诺保证使用ab47355于以下的经测试应用

“应用说明”部分 下显示的仅为推荐的起始稀释度;实际最佳的稀释度/浓度应由使用者检定。

应用	Ab评论	说明
Sandwich ELISA		Use at an assay dependent concentration.

图片



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.cn/abpromise> or contact our technical team.

### **Terms and conditions**

---

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