

Mouse VCAM1 ELISA Kit ab100750

5 References 2 图像

概述

产品名称	小鼠VCAM1 ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 1.5 pg/ml
范围	2.05 pg/ml - 500 pg/ml
回收率	96 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	97.36	84% - 107%
Serum	96.43	82% - 105%
Plasma	94.64	81% - 104%

实验步骤	Multiple steps standard assay
种属反应性	与反应: Mouse
产品概述	<p>Abcam's VCAM1 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of mouse VCAM1 in serum, plasma and cell culture supernatants.</p> <p>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.</p>
平台	Microplate

性能

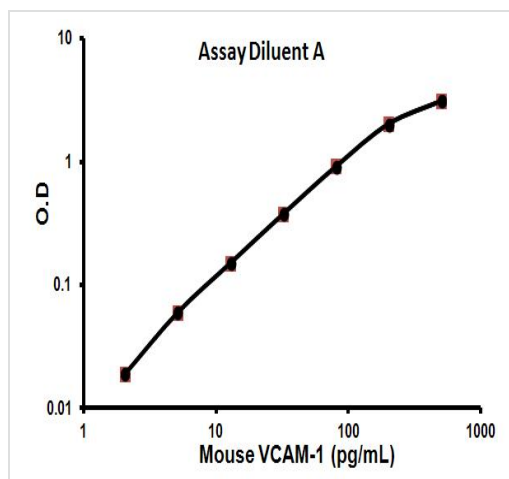
存放说明

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

组件	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

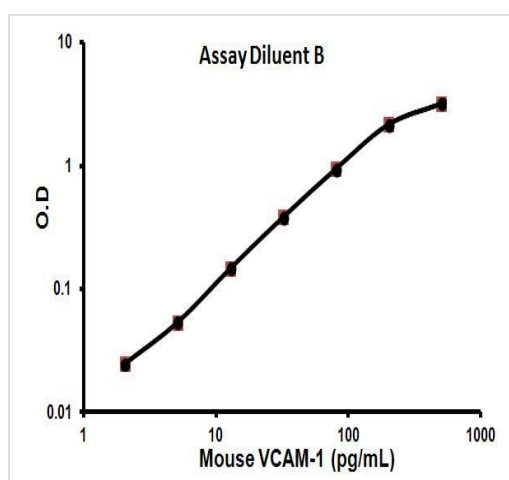
功能	Important in cell-cell recognition. Appears to function in leukocyte-endothelial cell adhesion. Interacts with the beta-1 integrin VLA4 on leukocytes, and mediates both adhesion and signal transduction. The VCAM1/VLA4 interaction may play a pathophysiologic role both in immune responses and in leukocyte emigration to sites of inflammation.
组织特异性	Expressed on inflamed vascular endothelium, as well as on macrophage-like and dendritic cell types in both normal and inflamed tissue.
序列相似性	Contains 7 Ig-like C2-type (immunoglobulin-like) domains.
结构域	Either the first or the fourth Ig-like C2-type domain is required for VLA4-dependent cell adhesion.
翻译后修饰	Sialoglycoprotein.
细胞定位	Membrane.

图片



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.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