

Human CD62E ELISA Set (without plates) ab47348

1 图像

概述

产品名称	人CD62E ELISA试剂盒(without plates)
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 0.5 ng/ml
范围	1 ng/ml - 32 ng/ml
检测时间	2h 15m
实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	Abcam's CD26E Human <i>in vitro</i> ELISA Set is designed for the quantitative measurement of CD26E in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.

The CD62E 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.

This ELISA set is intended to measure human CD62E 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.

Recognizes both natural and recombinant human CD62E.

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	标识符	10 x 96 tests	5 x 96 tests
CD62E Capture Antibody		2 x 0.5ml	1 x 0.5ml

组件	标识符	10 x 96 tests	5 x 96 tests
CD62E standard	Yellow	10 vials	5 vials
Chromogen TMB Substrate Solution		4 vials	2 vials
Detection biotinylated anti-CD62E antibody		2 x 0.55ml	1 x 0.55ml
Streptavidin-HRP		2 vials	1 vial

相关性 CD62E is expressed on cytokine induced endothelial cells and mediates their binding to leukocytes. The ligand recognized by CD62E is sialyl-lewis X (alpha(1->3)fucosylated derivatives of polylectosamine that are found at the nonreducing termini of glycolipids).

细胞定位 Membrane; single-pass type I membrane protein

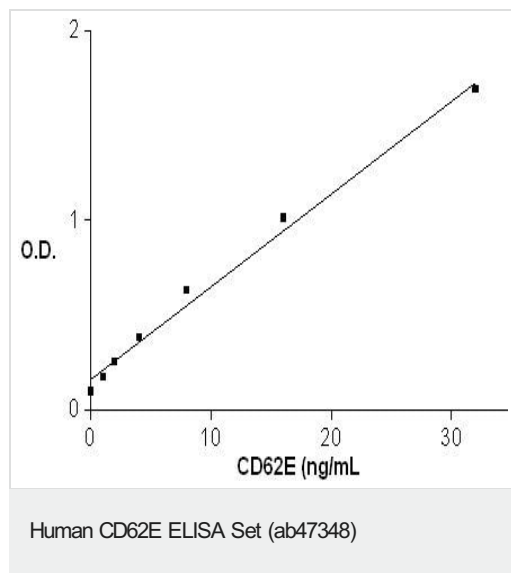
应用

The Abpromise guarantee **Abpromise™**承诺保证使用ab47348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



CD62E Standard curve ranging from 1 to 32 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