

### Human CD62E ELISA Kit (Selectin) ab100512

[1 References](#) [2 图像](#)

#### 概述

产品名称	人CD62E ELISA试剂盒(Selectin)
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 30 pg/ml
范围	24.69 pg/ml - 18000 pg/ml
回收率	94 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	94.39	83% - 103%
Serum	93.75	83% - 102%
Plasma	95.38	84% - 105%

**实验步骤** Multiple steps standard assay

**种属反应性** 与反应: Human

**产品概述** Abcam's CD62E (Selectin) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human CD62E in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human CD62E coated on a 96-well plate. Standards and samples are pipetted into the wells and CD62E present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human CD62E 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 CD62E bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

**说明** Optimization may be required with urine samples

**平台** Microplate

## 性能

### 存放说明

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

组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
900x HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated anti-Human CD62E	2 vials
CD62E Microplate (12 strips x 8 wells)	1 unit
Recombinant Human CD62E Standard (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

### 功能

Cell-surface glycoprotein having a role in immunoadhesion. Mediates in the adhesion of blood neutrophils in cytokine-activated endothelium through interaction with PSGL1/SELPLG. May have a role in capillary morphogenesis.

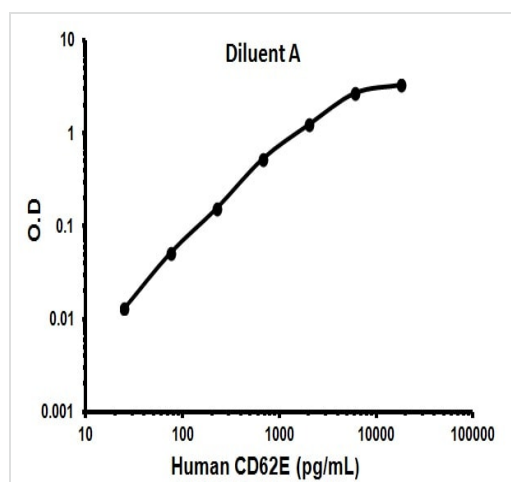
### 序列相似性

Belongs to the selectin/LECAM family.  
Contains 1 C-type lectin domain.  
Contains 1 EGF-like domain.  
Contains 6 Sushi (CCP/SCR) domains.

### 细胞定位

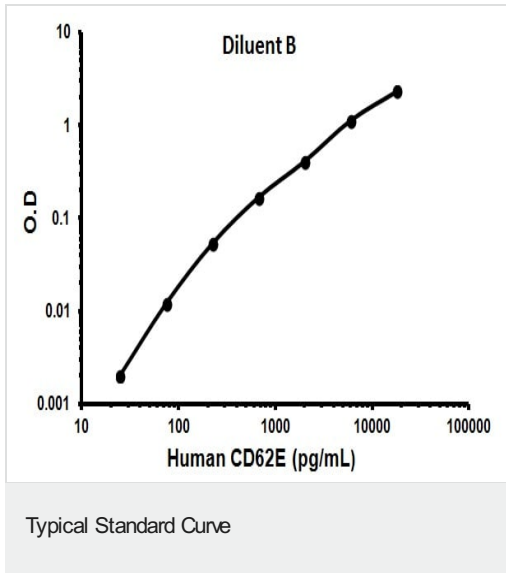
Membrane.

## 图片



Representative Standard Curve using ab100512.

Typical Standard Curve



Representative Standard Curve using ab100512.

**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