

Human CD62E ELISA Kit (Selectin) ab100512

1 References 画像数 2

製品の概要

製品名	Human CD62E ELISA Kit (Selectin)
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 30 pg/ml
検出範囲	24.69 pg/ml - 18000 pg/ml
添加回収試験	94 %

特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	94.39	83% - 103%
Serum	93.75	83% - 102%
Plasma	95.38	84% - 105%

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 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.

特記事項 Optimization may be required with urine samples

試験プラットフォーム Microplate

製品の特性

保存方法Store at -20°C. Please refer to protocols.

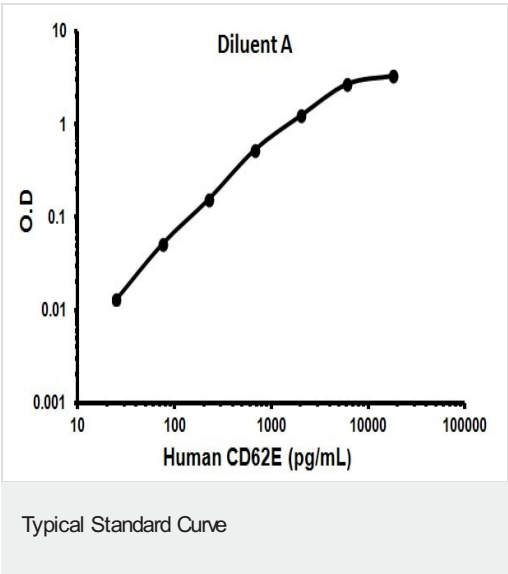
内容	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
900x HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated anti-Human CD62E	2 vials
CD62E Microplate (12 strips x 8 wells)	1 unit
Recombinant Human CD62E Standard (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB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

機能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.

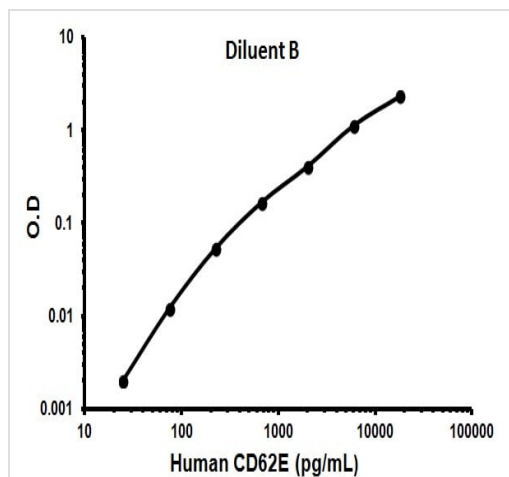
配列類似性Belongs to the selectin/LECAM family.
Contains 1 C-type lectin domain.
Contains 1 EGF-like domain.
Contains 6 Sushi (CCP/SCR) domains.

細胞内局在Membrane.

画像



Representative Standard Curve using ab100512.



Representative Standard Curve using ab100512.

Typical 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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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