

Rat CD62L ELISA Kit (Selectin L) ab100776

画像数 2

製品の概要

| | |
|---------|---|
| 製品名 | Rat CD62L ELISA Kit (Selectin L) |
| 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 | < 6 pg/ml |
| 検出範囲 | 8.23 pg/ml - 6000 pg/ml |
| 添加回収試験 | 96 % |

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 | 98.42 | 89% - 107% |
| Serum | 97.06 | 86% - 107% |
| Plasma | 94.33 | 85% - 103% |

ステップ Multiple steps standard assay

種交差性 交差種: Rat

製品の概要 Abcam's CD62L Rat ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Rat CD62L in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for CD62L coated on a 96-well plate. Standards and samples are pipetted into the wells and CD62L present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Rat CD62L 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 CD62L bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

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 |
| 500X HRP-Streptavidin Concentrate | 1 x 200µl |
| Assay Diluent A | 1 x 30ml |
| Biotinylated anti Rat CD62L | 2 vials |
| CD62L Microplate (12 x 8 wells) | 1 unit |
| Recombinant rat CD62L Standard (lyophilized) | 2 vials |
| Stop Solution | 1 x 8ml |
| TMB One-Step Substrate Reagent | 1 x 12ml |

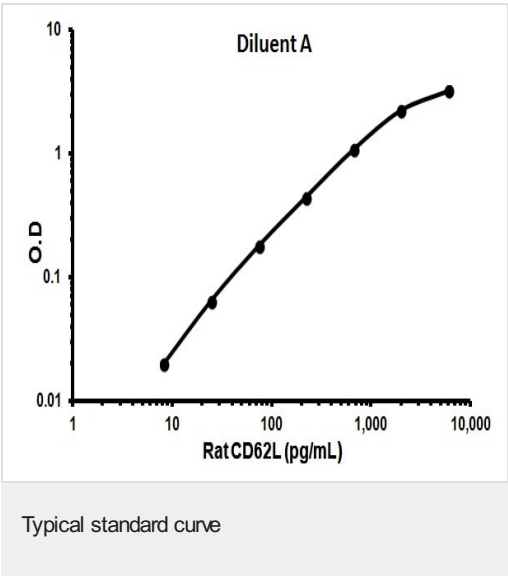
関連性

CD62L is a cell surface component that is a member of a family of adhesion/homing receptors. These play important roles in leukocyte-endothelial cell interactions. CD62L is composed of multiple domains: one homologous to lectins, one to epidermal growth factor, and two to the consensus repeat units found in C3/C4 binding proteins.

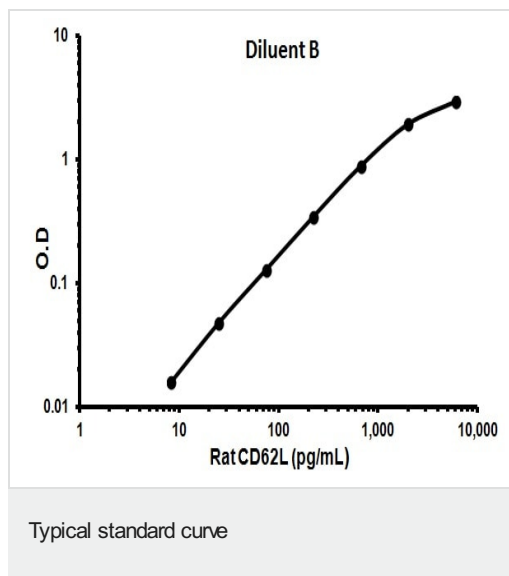
細胞内局在

Cell Membrane; single-pass type I membrane protein.

画像



Representative standard curve using ab100776



Representative standard curve using ab100776

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