

Rat CD62L ELISA Kit (Selectin L) ab100776

2 图像

概述

产品名称	大鼠CD62L ELISA试剂盒(Selectin L)
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 6 pg/ml
范围	8.23 pg/ml - 6000 pg/ml
回收率	96 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	98.42	89% - 107%
Serum	97.06	86% - 107%
Plasma	94.33	85% - 103%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Rat

产品概述 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.

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
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

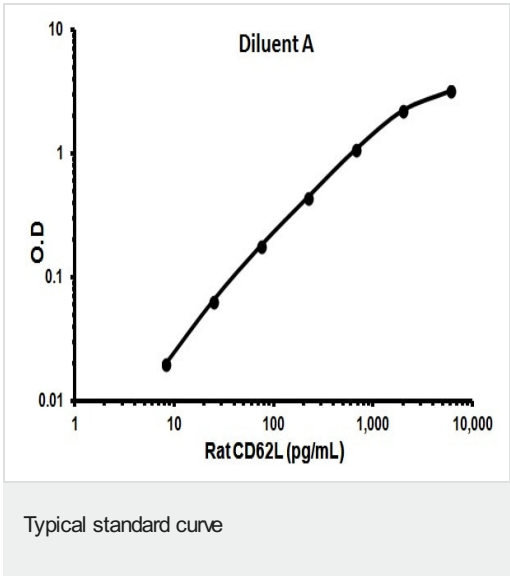
相关性

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.

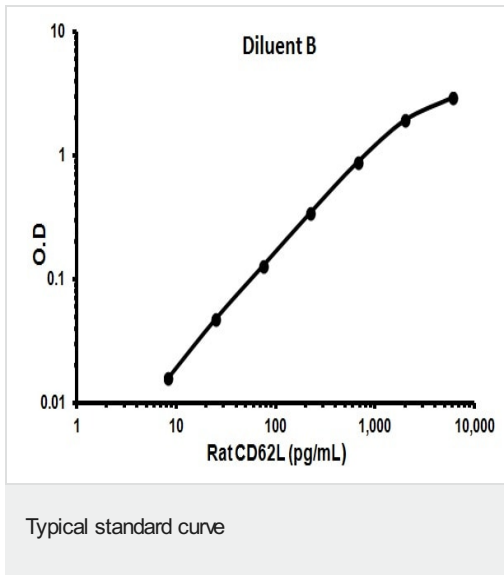
细胞定位

Cell Membrane; single-pass type I membrane protein.

图片



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.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