

Mouse CD62P ELISA Kit (P-Selectin) ab100737

画像数 2

製品の概要

製品名	Mouse CD62P ELISA Kit (P-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	< 8 pg/ml
検出範囲	8.23 pg/ml - 6000 pg/ml
添加回収試験	93 %

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	94.62	83% - 105%
Serum	92.38	80% - 102%
Plasma	93.57	81% - 104%

ステップ Multiple steps standard assay

種交差性 交差種: Mouse

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

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

Get higher sensitivity in only 90 minutes with Mouse P-Selectin ELISA Kit (CD62P) ([ab200014](#))

from our SimpleStep ELISA® range.

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
600X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated anti-mouse P-Selectin	2 vials
P-Selectin Microplate (12 x 8 wells)	1 unit
Recombinant Mouse P-Selectin Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

機能

Ca(2+)-dependent receptor for myeloid cells that binds to carbohydrates on neutrophils and monocytes. Mediates the interaction of activated endothelial cells or platelets with leukocytes. The ligand recognized is sialyl-Lewis X. Mediates rapid rolling of leukocyte rolling over vascular surfaces during the initial steps in inflammation through interaction with PSGL1.

組織特異性

Stored in the alpha-granules of platelets and Weibel-Palade bodies of endothelial cells. Upon cell activation by agonists, P-selectin is transported rapidly to the cell surface.

関連疾患

Defects in SELP may be a cause of susceptibility to ischemic stroke (ISCHSTR) [MIM:601367]; also known as cerebrovascular accident or cerebral infarction. A stroke is an acute neurologic event leading to death of neural tissue of the brain and resulting in loss of motor, sensory and/or cognitive function. Ischemic strokes, resulting from vascular occlusion, is considered to be a highly complex disease consisting of a group of heterogeneous disorders with multiple genetic and environmental risk factors.

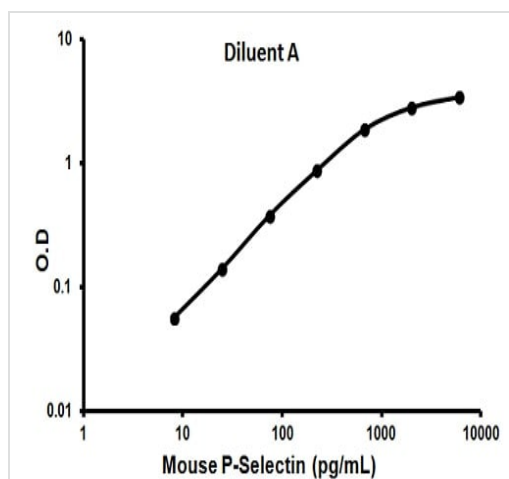
配列類似性

Belongs to the selectin/LECAM family.
Contains 1 C-type lectin domain.
Contains 1 EGF-like domain.
Contains 9 Sushi (CCP/SCR) domains.

細胞内局在

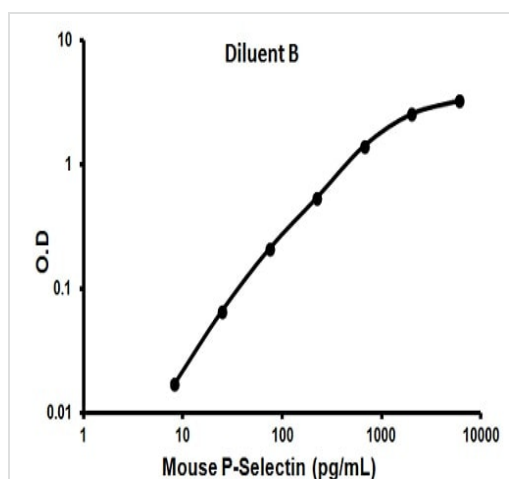
Membrane.

画像



Representative standard curve using ab100737

Typical standard curve



Representative standard curve using ab100737

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