

## 4 References 画像数 2

## 製品の概要

製品名	Human 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	< 20 pg/ml
検出範囲	0.041 ng/ml - 30 ng/ml
添加回収試験	72 %

### 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	72.51	63% - 85%
Serum	74.98	65% - 86%
Plasma	68.29	60% - 78%

ステップ	Multiple steps standard assay
種交差性	交差種: Human
製品の概要	<p>Abcam's CD62P (P-Selectin) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human CD62P in serum, plasma, and cell culture supernatants.</p> <p>This assay employs an antibody specific for Human CD62P coated on a 96-well plate. Standards and samples are pipetted into the wells and CD62P present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human CD62P 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 CD62P bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.</p>
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
800X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated anti-Human CD62P	2 vials
CD62P Microplate (12 x 8 wells)	1 unit
Recombinant Human CD62P 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.

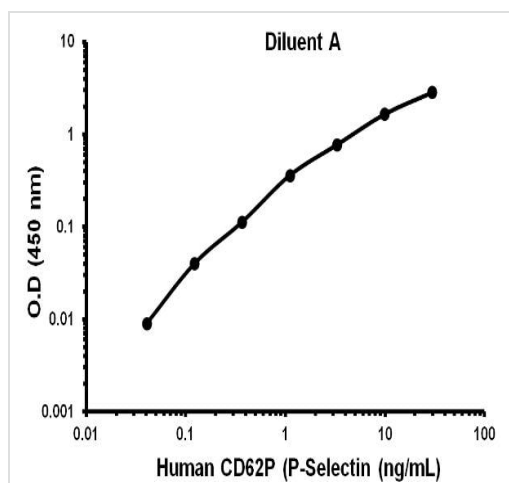
### 配列類似性

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

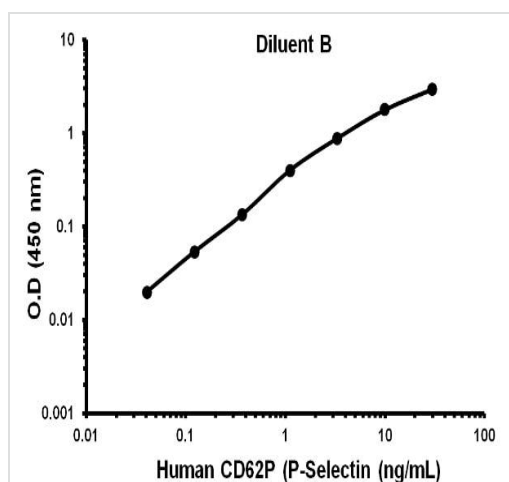
Membrane.

## 画像



Representative Standard Curve using ab100631.

Typical Standard Curve



Representative Standard Curve using ab100631.

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