

Product datasheet

Monkey CD62P ELISA Kit ab119554

画像数 1

製品の概要

製品名 Monkey CD62P ELISA Kit

検出方法 Colorimetric

再現性 Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
Overall	4			< 10%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Overall	4			< 10%

サンプルの種類 Cell culture supernatant, Serum

アッセイタイプ Sandwich (quantitative)

検出感度 1.3 U/ml

検出範囲 0.8 U/ml - 50 U/ml

添加回収試験

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
Serum	76	% - %

ステップ Multiple steps standard assay

種交差性 交差種: Monkey

製品の概要

Abcam's CD62P Monkey *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for accurate quantitative measurement of Monkey CD62P concentrations in cell culture supernatant, and serum (baboon).

CD62P specific antibody has been precoated onto 96-well plates. Standards and test samples are added to the wells and along with an HRP-conjugated CD62P detection antibody and the microplate is then incubated at room temperature. After the removal of unbound proteins by washing, TMB is used to visualize the HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changes into yellow after adding acidic stop solution. The

density of yellow coloration is directly proportional to the CD62P captured in the plate.

特記事項

Standard is recombinant protein produced in *E.coli*.

アプリケーション

適用あり: Sandwich ELISA

試験プラットフォーム

Microplate

製品の特性

保存方法

Store at +4°C. Please refer to protocols.

内容	1 x 96 tests
20X Assay Buffer Concentrate	1 x 5ml
20X Wash Buffer Concentrate	1 x 50ml
Adhesive Films	2 units
HRP-Conjugate anti-monkey CD62P monoclonal antibody	1 x 60µl
Microplate coated with monoclonal antibody to monkey CD62P (12 x 8 wells)	1 unit
Monkey CD62P Standard Lyophilized (100 U/mL upon reconstitution)	2 vials
Sample Diluent	1 x 12ml
Stop Solution (1M Phosphoric acid)	1 x 15ml
TMB Substrate Solution	1 x 15ml

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.

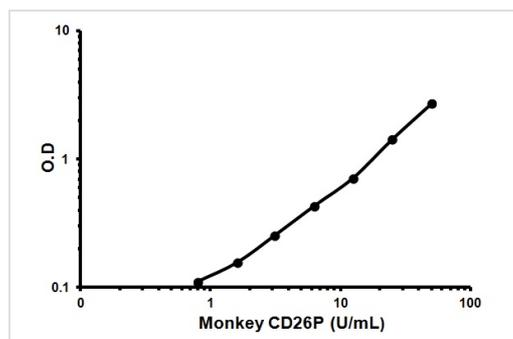
アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab119554** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

アプリケーション	Abreviews	特記事項
Sandwich ELISA		Use at an assay dependent concentration.

画像



Representative Standard Curve using ab119554.

Typical Standard Curve

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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