

Human CD62E ELISA Set (without plates) ab47348

画像数 1

製品の概要

製品名	Human CD62E ELISA Set (without plates)
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	< 0.5 ng/ml
検出範囲	1 ng/ml - 32 ng/ml
全工程の試験時間	2h 15m
ステップ	Multiple steps standard assay
種交差性	交差種: Human
製品の概要	<p>Abcam's CD26E Human <i>in vitro</i> ELISA Set is designed for the quantitative measurement of CD26E in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.</p> <p>The CD62E ELISA Set includes both the capture and the detection antibodies as well as Horse-Radish peroxidase conjugated Streptavidin (HRP-Strep) and TMB which are the reagents used in the colour development step.</p> <p>This ELISA set is intended to measure human CD62E in cell -culture supernatants, as well as in serum and plasma samples. The enclosed procedures provide general guidelines for ELISA preparation. This procedure may need some optimization depending on the plastic, matrix, standard range.</p> <p>Recognizes both natural and recombinant human CD62E.</p>
アプリケーション	適用あり: Sandwich ELISA
試験プラットフォーム	Reagents

製品の特性

保存方法 Store at +4°C. Please refer to protocols.

内容	ラベル	10 x 96 tests	5 x 96 tests
CD62E Capture Antibody		2 x 0.5ml	1 x 0.5ml

内容	ラベル	10 x 96 tests	5 x 96 tests
CD62E standard	Yellow	10 vials	5 vials
Chromogen TMB Substrate Solution		4 vials	2 vials
Detection biotinylated anti-CD62E antibody		2 x 0.55ml	1 x 0.55ml
Streptavidin-HRP		2 vials	1 vial

関連性 CD62E is expressed on cytokine induced endothelial cells and mediates their binding to leukocytes. The ligand recognized by CD62E is sialyl-lewis X (alpha(1->3)fucosylated derivatives of polylactosamine that are found at the nonreducing termini of glycolipids).

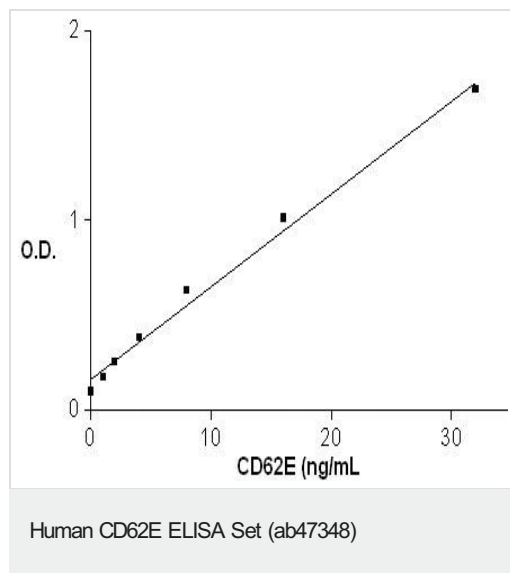
細胞内局在 Membrane; single-pass type I membrane protein

アプリケーション

The Abpromise guarantee Abpromise保証は、 次のテスト済みアプリケーションにおけるab47348の使用に適用されます
アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご検討ください。

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

画像



CD62E Standard curve ranging from 1 to 32 ng/mL

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