

Human CD62E ELISA Set (without plates) ab47348

1 Image

Overview

Product name	Human CD62E ELISA Set (without plates)
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 0.5 ng/ml
Range	1 ng/ml - 32 ng/ml
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Human
Product overview	<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>
Tested applications	Suitable for: Sandwich ELISA
Platform	Reagents

Properties

Storage instructionsStore at +4°C. Please refer to protocols.

Components	Identifier	10 x 96 tests	5 x 96 tests
CD62E Capture Antibody		2 x 0.5ml	1 x 0.5ml

Components	Identifier	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

Relevance

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).

Cellular localization

Membrane; single-pass type I membrane protein

Applications

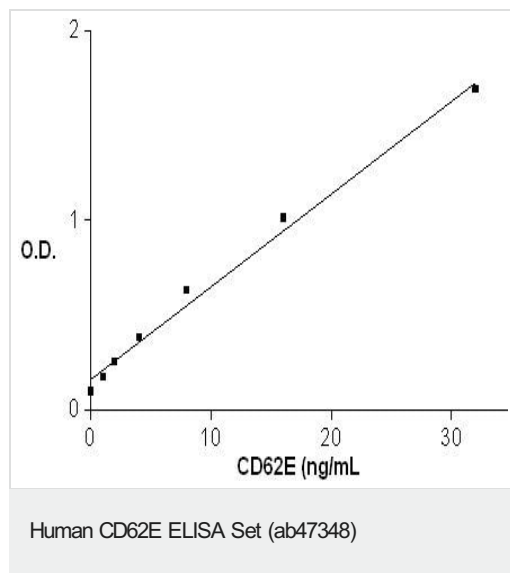
The Abpromise guarantee

Our **Abpromise guarantee** covers the use of ab47348 in the following tested applications.

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

Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

Images



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.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors