

Mouse L-Selectin ELISA Kit (CD62L) αb155448

1 References 1 Image

Overview

Product name	Mouse L-Selectin ELISA Kit (CD62L)				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 10%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 12%
Sample type	Cell culture supernatant, Serum, Plasma				
Assay type	Sandwich (quantitative)				
Sensitivity	< 35 pg/ml				
Range	40.96 pg/ml - 10000 pg/ml				
Recovery	Sample specific recovery				
	Sample type		Average %		Range
	Serum		103.1		95% - 111%
	Plasma		95.69		82% - 104%
	Cell culture media		109.4		103% - 116%
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Mouse				
Product overview	Abcam's L-Selectin (CD62L) Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay designed for the quantitative measurement of mouse L-Selectin in serum, plasma and cell culture supernatants.				

This assay employs an antibody specific for mouse L-Selectin (CD62L) coated on a 96-well plate.

Standards and samples are pipetted into the wells and L-Selectin present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse L-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 L-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 L-Selectin ELISA Kit (CD62L) ([ab199085](#)) from our SimpleStep ELISA® range.

Platform

Microplate

Properties**Storage instructions**

Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
700X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated anti-mouse L-Selectin	2 vials
L-Selectin Microplate (12 x 8 wells)	1 unit
Recombinant Mouse L-Selectin Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

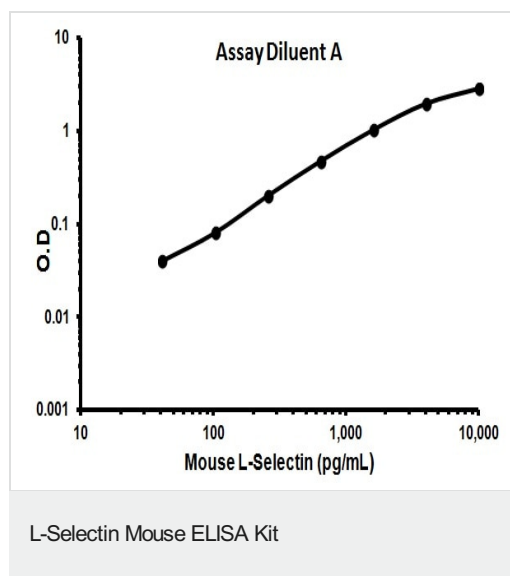
Relevance

CD62L is a cell surface component that is a member of a family of adhesion/homing receptors. These play important roles in leukocyte-endothelial cell interactions. CD62L is composed of multiple domains: one homologous to lectins, one to epidermal growth factor, and two to the consensus repeat units found in C3/C4 binding proteins.

Cellular localization

Cell Membrane; single-pass type I membrane protein.

Images



Representative standard curve using ab155448

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