

# Human CD62L ELISA Kit (Selectin L) ab45917

[1 References](#)   [1 Image](#)

Overview

Product name	Human CD62L ELISA Kit (Selectin L)				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Sample A	15	28ng/ml	1.3	4.6%
	Sample B	19	9ng/ml	0.5	5.5%
Sample type	Cell culture supernatant, Serum, Plasma				
Assay type	Sandwich (quantitative)				
Sensitivity	< 1 ng/ml				
Range	1.9 ng/ml - 60 ng/ml				
Assay time	1h 45m				
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Human				
Product overview	<p>Abcam's CD62L (Selectin L) Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of CD62L in Human serum, plasma, buffered solutions and cell culture medium.</p> <p>A monoclonal antibody specific for CD62L has been coated onto the wells of the microtiter strips provided. Samples, including standards of known CD62L concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for CD62L are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated and washed. A TMB substrate solution is added which acts on the bound enzyme to induce a colored reaction product. The intensity of this colored product is directly proportional to the concentration of CD62L present in the samples.</p> <p>This kit will recognize both endogenous and recombinant Human CD62L.</p>				
Platform	Microplate				

Properties

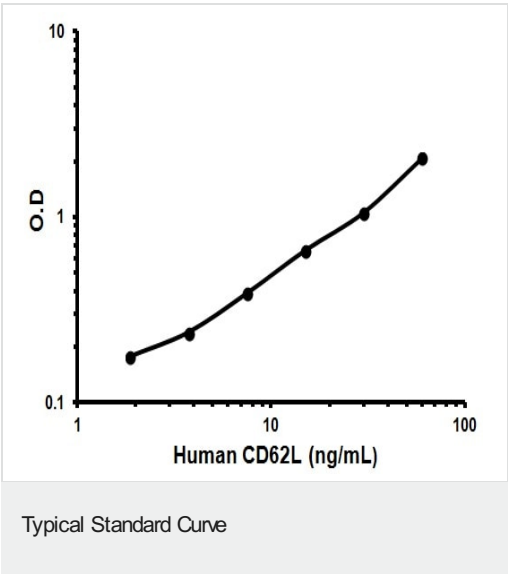
**Storage instructions** Store at +4°C. Please refer to protocols.

Components	Identifier	1 x 96 tests	2 x 96 tests
10X Standard Diluent Buffer	Black	1 x 15ml	1 x 25ml
200X Wash Buffer	White	1 x 10ml	2 x 10ml
Biotinylated Antibody Diluent	Red	1 x 7ml	1 x 13ml
Biotinylated anti-CD62L antibody	Red	1 x 0.4ml	2 x 0.4ml
CD62L Control	Silver	2 vials	4 vials
CD62L Microplate (12 x 8 well strips)		1 unit	2 units
CD62L Standard	Yellow	2 vials	4 vials
Chromogen TMB Substrate Solution		1 x 11ml	1 x 24ml
HRP Diluent	Red	1 x 11ml	1 x 23ml
Stop Reagent	Black	1 x 11ml	2 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl

**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 ab45917

**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