

# PRODUCT INFORMATION



## E-Selectin/CD62E Extracellular Domain (human, recombinant)

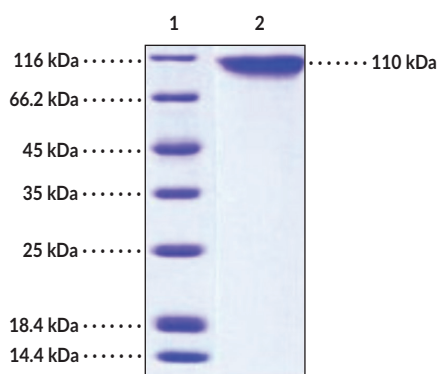
Item No. 35725

### Overview and Properties

**Synonyms:** CD62 Antigen-like Family Member E, ELAM1, Endothelial Leukocyte Adhesion Molecule 1, LECAM2, Leukocyte-endothelial Cell Adhesion Molecule 2  
**Source:** Recombinant human C-terminal His-tagged E-selectin expressed in HEK293 cells  
**Amino Acids:** 22-556  
**Uniprot No.:** P16581  
**Molecular Weight:** 60 kDa  
**Storage:** -80°C (as supplied)  
**Stability:** ≥1 year  
**Purity:** ≥97% estimated by SDS-PAGE  
**Supplied in:** Lyophilized from sterile PBS, pH 7.4  
**Endotoxin Testing:** <1.0 EU/μg, determined by the LAL endotoxin assay

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Image



Lane 1: MW Markers  
Lane 2: E-Selectin Extracellular Domain

**SDS-PAGE Analysis of E-Selectin Extracellular Domain.** This protein has a calculated molecular weight of 60 kDa. It has an apparent molecular weight of approximately 110 kDa by SDS-PAGE under reducing conditions due to glycosylation.

WARNING  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/08/2021

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

E-Selectin, also known as CD62E, is a glycoprotein encoded by the *SELE* gene in humans that has roles as a cellular adhesion molecule.<sup>1,2</sup> It is composed of an N-terminal C-type lectin-like domain, which recognizes carbohydrate ligands, an EGF-like domain, multiple short consensus repeats (SCRs), a transmembrane region, and a C-terminal cytoplasmic tail.<sup>2,3</sup> E-Selectin is constitutively expressed in bone marrow endothelial cells, where it has a role in maintaining the stem cell niche, and its expression is induced in vascular endothelial cells by inflammatory stimuli, such as TNF- $\alpha$ , IL-1 $\beta$ , or LPS.<sup>1,2</sup> It also exists as a soluble form that is shed from the cell surface *via* proteolytic cleavage.<sup>4</sup> E-selectin recognizes and binds to a variety of glycoproteins and glycolipids expressed by leukocytes, hematopoietic cells, or cancer cells, including L-selectin, sialyl Lewis X (sLe<sup>x</sup>), and E-selectin ligand-1 (ESL-1), as well as CD44 and CD43E.<sup>5</sup> Binding of E-selectin to its ligands facilitates the tethering and rolling of infiltrating cells on endothelial cells and promotes their extravasation into inflamed tissues.<sup>2</sup> *Sele*<sup>-/-</sup> mice exhibit reduced infiltration of metastatic cancer cells to the lung, and serum soluble E-selectin levels are elevated in patients with a variety of conditions, including atherosclerosis, type 1- and type 2 diabetes, sepsis, and multiple sclerosis.<sup>6-9</sup> Cayman's E-Selectin/CD62E Extracellular Domain (human, recombinant) protein consists of 546 amino acids, has a calculated molecular weight of 60 kDa, and a predicted N-terminus of Trp22 after signal peptide cleavage. By SDS-PAGE, under reducing conditions, the apparent molecular mass of the protein is 110 kDa due to glycosylation.

## References

1. Silva, M., Videira, P.A., and Sackstein, R. E-selectin ligands in the human mononuclear phagocyte system: Implications for infection, inflammation, and immunotherapy. *Front. Immunol.* **8**, 1878 (2018).
2. Tvaroška, I., Selvaraj, C., and Koča, J. Selectins-The two Dr. Jekyll and Mr. Hyde faces of adhesion molecules-A review. *Molecules* **25(12)**, 2835 (2020).
3. Aleisa, F.A., Sakashita, K., Lee, J.M., *et al.* Functional binding of E-selectin to its ligands is enhanced by structural features beyond its lectin domain. *J. Biol. Chem.* **295(11)**, 3719-3733 (2020).
4. DeLisser, H.M., Christofidou-Solomidou, M., Sun, J., *et al.* Loss of endothelial surface expression of E-selectin in a patient with recurrent infections. *Blood* **94(3)**, 884-894 (1999).
5. Barthel, S.R., Gavino, J.D., Descheny, L., *et al.* Targeting selectins and selectin ligands in inflammation and cancer. *Expert Opin. Ther. Targets* **11(11)**, 1473-1491 (2007).
6. Hiratsuka, S., Goel, S., Kamoun, W.S., *et al.* Endothelial focal adhesion kinase mediates cancer cell homing to discrete regions of the lungs via E-selectin up-regulation. *Proc. Natl. Acad. Sci. USA* **108(9)**, 3725-3730 (2011).
7. Paterson, A.D., Lopes-Virella, M.F., Waggott, D., *et al.* Genome-wide association identifies the ABO blood group as a major locus associated with serum levels of soluble E-selectin. *Arterioscler. Thromb. Vasc. Biol.* **29(11)**, 1958-1967 (2009).
8. Cummings, C.J., Sessler, C.N., Beall, L.D., *et al.* Soluble E-selectin levels in sepsis and critical illness. Correlation with infection and hemodynamic dysfunction. *Am. J. Respir. Crit. Care Med.* **156(2 Pt 1)**, 431-437 (1997).
9. Tsukada, N., Miyagi, K., Matsuda, M., *et al.* Soluble E-selectin in the serum and cerebrospinal fluid of patients with multiple sclerosis and human T-lymphotropic virus type 1-associated myelopathy. *Neurology* **45(10)**, 1914-1918 (1995).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM