

PRODUCT INFORMATION



ICAM-1/CD54 Extracellular Domain (human, recombinant)

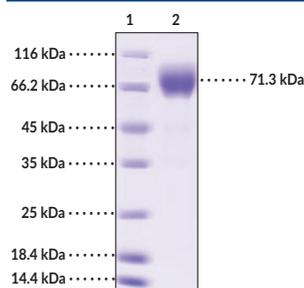
Item No. 44187

Overview and Properties

Synonyms: ICAM-1, Intercellular Adhesion Molecule-1
Source: Recombinant human C-terminal DDDDK-tagged ICAM-1 extracellular domain expressed in HEK293 cells
Amino Acids: 1-480
Uniprot No.: P05362
Molecular Weight: 50.2 kDa
Storage: -80°C (as supplied)
Stability: ≥1 year
Purity: ≥95% 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
Bioactivity: See figures for details

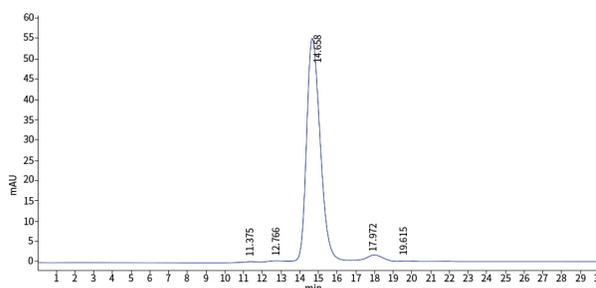
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: MW Markers
Lane 2: ICAM-1/CD54 Extracellular Domain

SDS-PAGE Analysis of ICAM-1/CD54 Extracellular Domain. This protein has a calculated molecular weight of 50.2 kDa. It has an apparent molecular weight of approximately 71.3 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, 12/23/2025

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

Intercellular adhesion molecule-1 (ICAM-1), also known as CD54, is a cell surface glycoprotein and member of the Ig superfamily.¹ It is composed of five extracellular Ig domains, which can vary in number through alternative splicing, a transmembrane domain, and a short C-terminal cytoplasmic domain.^{2,3} ICAM-1 is basally expressed at low levels but cytokines, such as NF- κ B, TNF- α , and IL-1 α , increase ICAM-1 levels in endothelial, epithelial, and immune cells.⁴ It binds integrins, such as lymphocyte function-associated antigen-1 (LFA-1) and complement 3 receptor (C3R), and directs lymphocytes or other immune cells to locally inflamed areas, among other roles in forming epithelial and endothelial barriers, regulating cell migration, and immune signaling.^{1,2,5} ICAM-1 can be cleaved from the cell membrane by cathepsin G or neutrophil elastase and promote or antagonize inflammation in a context-dependent manner and induce wound healing.^{1,3} Blood levels of soluble ICAM-1 are increased in patients with cancer, and serum levels of soluble ICAM-1 are increased in patients with acute ischemic stroke.^{6,7} Cayman's ICAM-1/CD54 Extracellular Domain (human, recombinant) protein consists of 459 amino acids and has a calculated molecular weight of 50.2 kDa. By SDS-PAGE, under reducing conditions, the apparent molecular mass of the protein is 71.3 kDa due to glycosylation.

References

1. Bui, T.M., Wiesolek, H.L. and Sumagin, R. ICAM-1: A master regulator of cellular responses in inflammation, injury resolution, and tumorigenesis. *J. Leukoc. Biol.* **108(3)**, 787-799 (2020).
2. Yang, Y., Jun, C.-D., Liu, J.-H., *et al.* Structural basis for dimerization of ICAM-1 on the cell surface. *Mol. Cell* **14(2)**, 269-276 (2004).
3. Robledo, O., Papaioannou, A., Ochietti, B., *et al.* ICAM-1 isoforms: Specific activity and sensitivity to cleavage by leukocyte elastase and cathepsin G. *Eur. J. Immunol.* **33(5)**, 1351-1360 (2003).
4. Yamada, Y., Tashiro, E., Taketani, S., *et al.* Mycotrienin II, a translation inhibitor that prevents ICAM-1 expression induced by pro-inflammatory cytokines. *J. Antibiot. (Tokyo)* **64(5)**, 361-366 (2011).
5. Yang, S.F., Chen, M.K., Hsieh, Y.S., *et al.* Prostaglandin E₂/EP1 signaling pathway enhances intercellular adhesion molecule 1 (ICAM-1) expression and cell motility in oral cancer cells. *J. Biol. Chem.* **285(39)**, 29808-29816 (2010).
6. Banks, R.E., Gearing, A.J.H., Hemingway, I.K., *et al.* Circulating intercellular adhesion molecule-1 (ICAM-1), E-selectin and vascular cell adhesion molecule-1 (VCAM-1) in human malignancies. *Br. J. Cancer* **68(1)**, 122-124 (1993).
7. Shyu, K.G., Chang, H. and Lin, C.C. Serum levels of intracellular adhesion molecule-1 and E-selectin in patients with acute ischaemic stroke. *J. Neurol.* **244(2)**, 90-93 (1997).