

PRODUCT INFORMATION



HVEM/CD270 Extracellular Domain (human, recombinant)

Item No. 31840

Overview and Properties

Synonyms: ATAR, Herpes Virus Entry Mediator, HVEA, LIGHTR, TNFRSF14, TR2, Tumor Necrosis Factor Receptor Superfamily Member 14

Source: Active recombinant human C-terminal His-tagged HVEM expressed in HEK293 cells

Amino Acids: 37-202

Uniprot No.: Q92956

Molecular Weight: 19 kDa

Storage: -80°C (as supplied)

Stability: ≥1 year

Purity: ≥90% estimated by SDS-PAGE

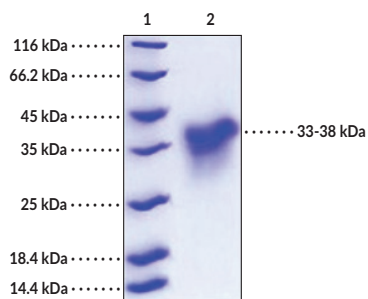
Supplied in: Lyophilized from sterile PBS, pH 7.4

Endotoxin Testing: <1.0 EU/μg, determined by the LAL endotoxin assay

Bioactivity: See figures for details

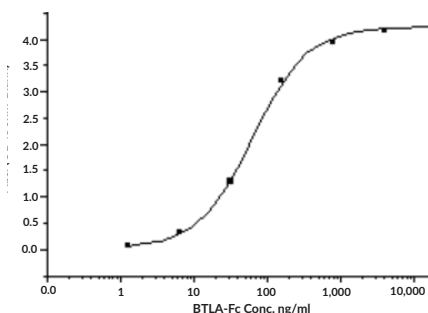
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: MW Markers
Lane 2: HVEM/CD270 Extracellular Domain

SDS-PAGE Analysis of HVEM/CD270 Extracellular Domain.
This protein has a calculated molecular weight of 19 kDa. It has an apparent molecular weight of approximately 33-38 kDa by SDS-PAGE under reducing conditions due to glycosylation.



HVEM/CD270 Extracellular Domain Binding in a Binding Assay.
Immobilized human HVEM/CD270 Extracellular Domain at 10 μg/ml (100 μl/well) can bind human BTLA-Fc with a linear range of 1.28-20 μg/ml.

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, 04/08/2021

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

Herpes virus entry mediator (HVEM), also known as CD270, is a type I transmembrane glycoprotein encoded by *TNFRSF14* in humans.^{1,2} It is composed of an extracellular domain containing four cysteine-rich repeats (CRDs), a transmembrane domain, and a cytoplasmic domain, which facilitates intracellular signaling. HVEM is expressed on T and B cells, other hematopoietic cells, and mucosal epithelium.² It binds to B and T lymphocyte attenuator (BTLA), CD160, and LIGHT (Item No. 31841), also known as CD258, and induces co-activation or co-inhibition of T cells in a ligand-dependent manner via signaling through TNF receptor-associated factors (TRAFs) or inhibitory tyrosine-based motifs (ITIMs), respectively.^{2,3} *Tnfrsf14*^{-/-} mice exhibit decreased susceptibility to bacterial infection, increased pro-inflammatory cytokine secretion and mortality in a model of concanavalin A-induced hepatitis, and increased susceptibility to the induction of experimental autoimmune encephalomyelitis (EAE), indicating roles for HVEM in both infection immunity and autoimmunity.⁴ Cayman's HVEM/CD270 Extracellular Domain (human, recombinant) protein can be used for ELISA and binding assay applications. This protein consists of 177 amino acids and has a calculated molecular weight of 19 kDa. By SDS-PAGE, under reducing conditions, the apparent molecular mass of the protein is approximately 33 to 38 kDa due to glycosylation.

References

1. Connolly, S.A., Landsburg, D.J., Carfi, A., *et al.* Structure-based analysis of the herpes simplex virus glycoprotein D binding site present on herpesvirus entry mediator HveA (HVEM). *J. Virol.* **76(21)**, 10894-10904 (2002).
2. Ware, C.F. Targeting the LIGHT-HVEM pathway. *Therapeutic targets of the TNF superfamily*. Grewal, I.S., editor, 1st ed., Springer (2009).
3. Shui, J.-W. and Kronenberg, M. HVEM: An unusual TNF receptor family member important for mucosal innate immune responses to microbes. *Gut Microbes* **4(2)**, 146-151 (2013).
4. Shui, J.-W., Steinberg, M.W., and Kronenberg, M. Regulation of inflammation, autoimmunity, and infection immunity by HVEM-BTLA signaling. *J. Leukoc. Biol.* **89(4)**, 517-523 (2011).

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