

PRODUCT INFORMATION

4-1BB Ligand Extracellular Domain (human, recombinant)

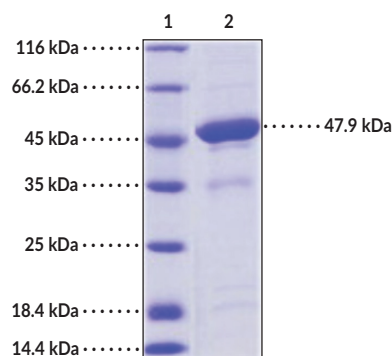
Item No. 32067

Overview and Properties

Synonyms:	4-1BBL, CD137L, CD137 Ligand, TNFSF9
Source:	Active recombinant N-terminal human IgG1 Fc-tagged 4-1BB ligand expressed in HEK293 cells
Amino Acids:	71-254
Uniprot No.:	P41273
Molecular Weight:	47.9 kDa
Storage:	-80°C (as supplied)
Stability:	≥1 year
Purity:	≥85% 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

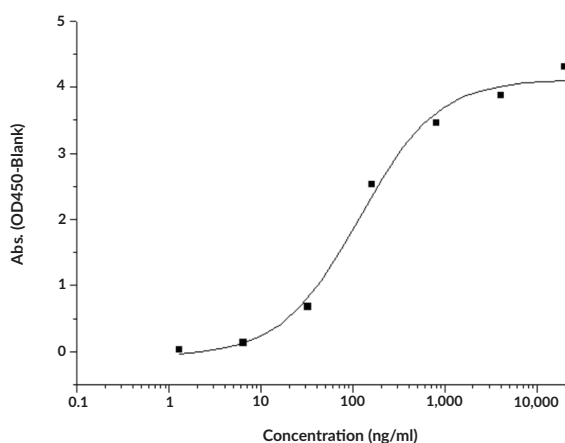
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: MW Markers
Lane 2: 4-1BB Ligand Extracellular Domain

SDS-PAGE Analysis of 4-1BB Ligand Extracellular Domain.



4-1BB Ligand Extracellular Domain Binding in a Functional ELISA.
Immobilized 4-1BB at 10 μg/ml can bind 4-1BB Ligand Extracellular Domain, the EC₅₀ value for this effect is 7-40 ng/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, 06/22/2020

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

TNF ligand superfamily member 9 (4-1BB ligand) is a type II transmembrane glycoprotein and a member of the TNF superfamily.^{1,2} It is composed of an N-terminal cytoplasmic region, a transmembrane domain, and a C-terminal extracellular tail region that contains a TNF homology domain.^{2,3} 4-1BB ligand production is induced in antigen-presenting cells (APCs), such as dendritic cells, B cells, and macrophages, by CD40 and toll-like receptor (TLR) activation.² It forms homotrimers and binds to its receptor, 4-1BB, on antigen-primed T cells, which induces NF- κ B and MAPK signaling through TNF receptor-associated factors (TRAFs) and, when the T cell is co-stimulated by other factors, induces PI3K-Akt signaling.^{1,2} This action promotes the proliferation of T cells, production of cytokines, and cell survival.² Soluble forms of 4-1BB ligand, consisting of only the extracellular domain, also form trimers that bind to its receptor but do not induce signaling.⁴ Mice deficient for 4-1BB ligand, *Tnfsf9*, have impairments in the formation of CD8⁺ T cell memory cells and in clearing pathogens.^{2,5} *Tnfsf9* knockout mice also develop B cell lymphomas before 12 months of age.⁶ 4-1BB ligand antibodies induce T regulatory cell expansion and prevent allogeneic graft rejection of pancreatic islets in mice.⁷ Intratumoral adenovirus-mediated gene transfer of *Tnfsf9* induces tumor regression in a mouse syngeneic model of breast cancer.⁶ The protein levels of 4-1BB ligand are elevated in more than 40% of patients with acute myeloid leukemia (AML) or non-Hodgkin's lymphoma but the expression of *TNFSF9* is either elevated or reduced in cancer patients depending on the type of cancer. Cayman's 4-1BB Ligand Extracellular Domain (human, recombinant) protein can be used for ELISA and binding assay applications. This protein is a disulfide-linked homodimer. The reduced monomer, comprised of 4-1BB ligand (amino acids 71-254) fused to human IgG1 Fc at the N-terminus, consists of 444 amino acids and has a calculated molecular weight of 47.9 kDa.

References

1. Barsoumian, H.B., Batra, L., Shrestha, P., *et al.* A novel form of 4-1BBL prevents cancer development via nonspecific activation of CD4⁺ T and natural killer cells. *Cancer Res.* **79**(4), 783-794 (2019).
2. So, T. and Ishii, N. The TNF-TNFR family of co-signal molecules. *Co-signal molecules in T cell activation.* Azuma, M. and Yagita, H., editors, 1st edition, *Springer Singapore* (2019).
3. Bitra, A., Doukov, T., Croft, M., *et al.* Crystal structures of the human 4-1BB receptor bound to its ligand 4-1BBL reveal covalent receptor dimerization as a potential signaling amplifier. *J. Biol. Chem.* **293**(26), 9958-9969 (2018).
4. Wyzgol, A., Müller, N., Fick, A., *et al.* Trimer stabilization, oligomerization, and antibody-mediated cell surface immobilization improve the activity of soluble trimers of CD27L, CD40L, 41BBL, and glucocorticoid-induced TNF receptor ligand. *J. Immunol.* **183**(3), 1851-1861 (2009).
5. Wang, C., Lin, G.H.Y., McPherson, A.J., *et al.* Immune regulation by 4-1BB and 4-1BBL: Complexities and challenges. *Immunol. Rev.* **229**(1), 192-215 (2009).
6. Vinay, D.S. and Kwon, B.S. Immunotherapy of cancer with 4-1BB. *Mol. Cancer Ther.* **11**(5), 1062-1070 (2012).
7. Elpek, K.G., Yolcu, E.S., Franke, D.D.H., *et al.* Ex vivo expansion of CD4⁺CD25⁺FoxP3⁺ T regulatory cells based on synergy between IL-2 and 4-1BB signaling. *J. Immunol.* **179**(11), 7295-7304 (2007).

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