

PRODUCT INFORMATION



B7-1/CD80 Long Isoform Extracellular Domain (human, recombinant; His- and Fc-tagged)

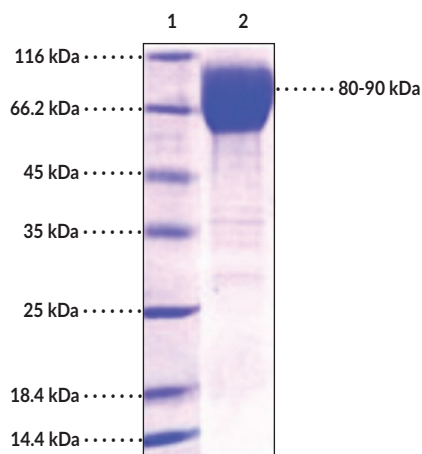
Item No. 31821

Overview and Properties

Synonyms: B7.1, BB1, CD28LG, CD28LG1, LAB7, T-lymphocyte Activation Antigen CD80
Source: Active recombinant C-terminal human IgG1 Fc-His-tagged CD80 expressed in HEK293 cells
Amino Acids: 35-242
Uniprot No.: P33681
Molecular Weight: 51.7 kDa
Storage: -20°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
Protein Concentration: *batch specific* mg/ml

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

Image



Lane 1: MW Markers

Lane 2: B7-1/CD80 Long Isoform Extracellular Domain (human, recombinant; His- and Fc-tagged)

SDS-PAGE Analysis of B7-1/CD80 Long Isoform Extracellular Domain (human, recombinant; His- and Fc-tagged). This protein has a calculated molecular weight of 51.7 kDa. It has an apparent molecular weight of approximately 80-90 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, 08/25/2022

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

CD80, also known as B7-1, is a glycoprotein and member of the CD28/B7 family of co-stimulatory receptors that promotes T cell activation.^{1,2} Alternative splicing of CD80 produces one full-length long isoform, CD80 long, and two short isoform, s1CD80 and s2CD80.² CD80 long exists as a membrane-bound dimer and is composed of an extracellular immunoglobulin variable (IgV) domain that interacts with the co-stimulatory molecule CD28 or the inhibitory molecule CTLA-4, as well as an immunoglobulin constant (IgC) domain and a cytoplasmic tail that are both required for T cell co-stimulation.¹ s1CD80 and s2CD80 lack the transmembrane domain or the transmembrane domain and the IgC domain, respectively, and are found in the serum.² CD80 is transiently expressed on the surface of antigen-presenting cells (APCs) and is upregulated following ligation of the co-stimulatory molecule CD86, as well as by pro-inflammatory stimuli, such as LPS, and downregulated by the anti-inflammatory cytokine IL-10.³⁻⁵ CD80 has two ligands, CD28 and CTLA-4, that compete for binding and are each expressed on naïve T cells but have opposing functions on T cell activation.⁶ CD80 binding to CD28 promotes T cell activation, survival, and cytokine production, whereas CD80 binding to CTLA-4 inhibits T cell activation and promotes T cell anergy. Neutralization of CD80 with a monoclonal antibody reduces the severity of synovitis and bone erosion, as well as CD4⁺ T cell infiltration, in inflamed joints in a mouse model of arthritis induced by complete Freund's adjuvant (CFA) and BSA.⁷ Decreased CD80 tumor levels have been identified in patients with renal cell carcinoma.⁸ CD80 SNPs have been found in individuals with gastric cancer.⁹ Cayman's B7-1/CD80 Long Isoform Extracellular Domain (human, recombinant; His- and FC-tagged) protein can be used for ELISA applications. This protein is a disulfide-linked homodimer. The reduced monomer, comprised of B7-1/CD80 (amino acids 35-242) fused to His-tagged human IgG1 Fc at its C-terminus, consists of 455 amino acids, a calculated molecular weight of 51.7 kDa, and a predicted N-terminus of Val35 after signal peptide cleavage. As a result of glycosylation, the monomer migrates at approximately 80-90 kDa by SDS-PAGE under reducing conditions.

References

1. Girard, T., Gaucher, D., El-Far, M., *et al.* CD80 and CD86 IgC domains are important for quaternary structure, receptor binding and co-signaling function. *Immunol. Lett.* **161(1)**, 65-75 (2014).
2. Kakoulidou, M., Giscombe, R., Zhao, X., *et al.* Human soluble CD80 is generated by alternative splicing, and recombinant soluble CD80 binds to CD28 and CD152 influencing T-cell activation. *Scand. J. Immunol.* **66(5)**, 529-537 (2007).
3. Mir, M.A. Costimulation in lymphomas and cancers. *Developing Costimulatory Molecules for Immunotherapy of Diseases*, Academic Press, Cambridge, MA (2015).
4. Sahoo, N.C., Rao, K.V.S., and Natarajan, K. CD80 expression is induced on activated B cells following stimulation by CD86. *Scand. J. Immunol.* **55(6)**, 577-584 (2002).
5. Zhou, F., Ciric, B., Li, H., *et al.* IL-10 deficiency blocks the ability of LPS to regulate expression of tolerance-related molecules on dendritic cells. *Eur. J. Immunol.* **42(6)**, 1449-1458 (2012).
6. Buchbinder, E.I. and Desai, A. CTLA-4 and PD-1 pathways: Similarities, differences, and implications of their inhibition. *Am. J. Clin. Oncol.* **39(1)**, 98-106 (2016).
7. Odobasic, D., Leech, M.T., Xue, J.R., *et al.* Distinct *in vivo* roles of CD80 and CD86 in the effector T-cell responses inducing antigen-induced arthritis. *Immunol. Invest.* **124(4)**, 503-513 (2008).
8. Flörcken, A., Johannsen, M., Nguyen-Hoai, T., *et al.* Immunomodulatory molecules in renal cell cancer: CD80 and CD86 are expressed on tumor cells. *Int. J. Clin. Exp. Pathol.* **10(2)**, 1443-1454 (2017).
9. Wu, R., Li, F., Zhu, J., *et al.* A functional variant at miR-132-3p, miR-212-3p, and miR-361-5p binding site in CD80 gene alters susceptibility to gastric cancer in a Chinese Han population. *Med. Oncol.* **31(8)**, 60 (2014).

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