

# PRODUCT INFORMATION



## ICOS/CD278 Extracellular Domain (human, recombinant)

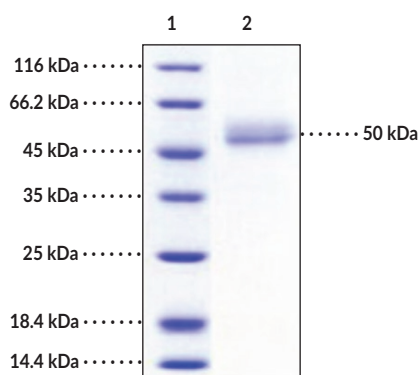
Item No. 32084

### Overview and Properties

<b>Synonyms:</b>	Activation-inducible Lymphocyte Immunomediatory Molecule, AILIM, CVID1, Inducible T Cell Costimulator
<b>Source:</b>	Active recombinant C-terminal human IgG1 Fc-His-tagged ICOS expressed in HEK293 cells
<b>Amino Acids:</b>	21-141
<b>Uniprot No.:</b>	Q9Y6W8
<b>Molecular Weight:</b>	41.6 kDa
<b>Storage:</b>	-80°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Purity:</b>	≥90% estimated by SDS-PAGE
<b>Supplied in:</b>	Lyophilized from sterile PBS, pH 7.4
<b>Endotoxin Testing:</b>	<1.0 EU/μg, determined by the LAL endotoxin assay
<b>Bioactivity:</b>	See figures for details

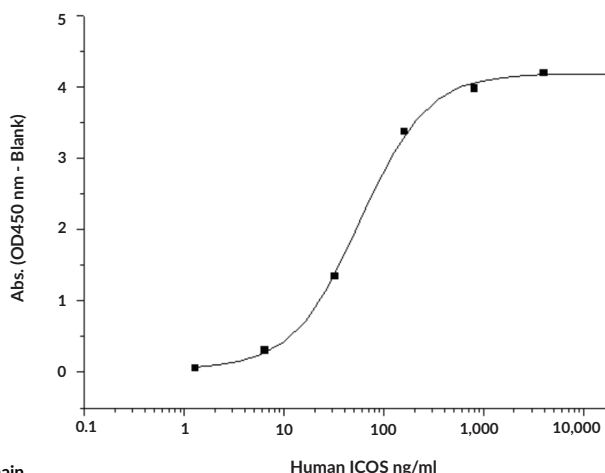
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Images



Lane 1: MW Markers  
Lane 2: ICOS Extracellular Domain

**SDS-PAGE Analysis of ICOS Extracellular Domain.**  
This protein has a calculated molecular weight of 41.6 kDa. It has an apparent molecular weight of approximately 50 kDa by SDS-PAGE under reducing conditions due to glycosylation.



**ICOS Bioactivity in functional ELISA.** Immobilized human B7-H2 at 1 μg/ml (100 μl/well) can bind human ICOS with a linear range of 1.6-200 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, 06/29/2020

**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

---

Inducible T cell co-stimulator (ICOS) is a glycoprotein and member of the CD28/B7 family of co-stimulatory receptors that promotes T cell activation.<sup>1</sup> It exists as a membrane-bound disulfide-linked homodimer and contains a leader peptide, an extracellular domain that interacts with ICOS ligand (ICOSL), and a cytoplasmic tail that mediates association with PI3K and is critical for T cell activation.<sup>1,2</sup> ICOS is expressed on activated CD4<sup>+</sup> and CD8<sup>+</sup> T cells following T cell receptor crosslinking and/or co-stimulation by CD28.<sup>1</sup> ICOS co-stimulation of T cells induces production of IL-4, IL-5, IL-6, IFN- $\gamma$  (Item Nos. 32008 | 32078), TNF- $\alpha$  (Item Nos. 32020 | 32069), and GM-CSF (Item Nos. 32044 | 32079), but not IL-2, and superinduces production of IL-10, as well as prevents T cell apoptosis.<sup>2</sup> Human melanoma cells express ICOSL and promote the expansion of ICOS<sup>+</sup> IL-10-secreting regulatory T cells, which is positively correlated with unfavorable clinical outcomes.<sup>3</sup> Administration of an anti-ICOS monoclonal antibody ameliorates mucosal inflammation and damage in a mouse model of T helper cell-induced chronic colitis.<sup>4</sup> Cayman's ICOS/CD278 Extracellular Domain (human, recombinant) protein can be used for ELISA. This protein is a disulfide-linked homodimer. The reduced monomer, comprised of ICOS (amino acids 21-141) fused to His-tagged human IgG1 Fc at its C-terminus, consists of 386 amino acids, has a calculated molecular weight of 41.6 kDa, and a predicted N-terminus of Glu21 after signal peptide cleavage. As a result of glycosylation, the monomer migrates at approximately 50 kDa by SDS-PAGE under reducing conditions.

## References

---

1. Wikenheiser, D.J. and Stumhofer, J.S. ICOS Co-Stimulation: Friend or Foe? *Front. Immunol.* **7**, 304 (2016).
2. Beier, K.C., Hutloff, A., Dittrich, A.M., *et al.* Induction, binding specificity and function of human ICOS. *Eur. J. Immunol.* **30(12)**, 3707-3717 (2000).
3. Amatore, F., Gorvel, L., and Olive, D. Inducible co-stimulator (ICOS) as a potential therapeutic target for anti-cancer therapy. *Expert Opin. Ther. Targets* **22(4)**, 343-351 (2018).
4. Kanai, T., Totsuka, T., Tezuka, K., *et al.* ICOS costimulation in inflammatory bowel disease. *J. Gastroenterol.* **37(Suppl 14)**, 78-81 (2002).