

## Data Sheet

### ICOS (CD278), Fc Fusion

Human, recombinant, Fc fusion protein

**Catalog#:** 71179

**Lot#:** 140722-D      **Conc.:** 1.54 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol

**Stability:** >6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Protein may be diluted to  $\geq 100 \mu\text{g/ml}$  in PBS + glycerol and stored at  $-80^{\circ}\text{C}$ .

**Endotoxin Level:** <1 EU/ $\mu\text{g}$

**References:**

1. Hutloff, A. *et al.*, *Nature*. 1999; **397**: 263-266.
2. Dong, C. *et al.*, *Nature*. 2001; **409**: 97-101.

**Description:**

Human ICOS, Fc fusion protein, also known as inducible T-cell costimulator and CD antigen 278, activation-inducible lymphocyte immunomediatory molecule, AILIM, CD278, and CVID1. GenBank Accession No. NM\_012092, a.a. 21-140 expressed in a HEK293 cell expression system. MW = 40.2 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

**Application:**

Useful for studying protein binding and for screening small molecules and antibodies.

### Quality Assurance

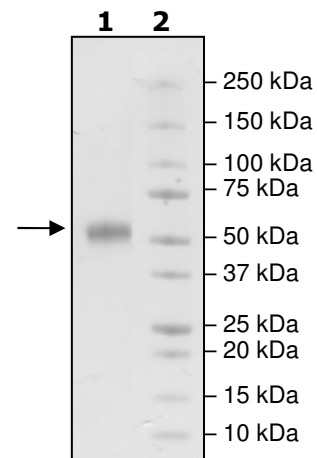
**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
3  $\mu\text{g}$  ICOS (CD278)

**Lane 2:**  
Protein Marker

**MW:** 40.2 kDa + glycans

**Purity:**  $\geq 90\%$



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6044 Cornerstone Court W, Ste E  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## ***Sequence***

**EINGSANYEMFIFHNGGVQILCKYPDIVQQFKMQLLKGGQILCDLTKTKGSGNTVSIKS  
LKFCHSQLSNNSVSFFLYNLDHSHANYFCNLSIFDPPPFKVTLTGGYLHIYESQLCCQ  
LKI**EGRMD**PKSCDKTHTCPPCPAPELLGGPSVFLFPPKPKDTLMISRTPEVTCVVDVS  
HEDPEVKFNWYVDGVEVHNAKTKPREEQYNSTYRVVSVLTVLHQDWLNGKEYKCKVS  
NKALPAIEKTISKAKGQPREPQVYTLPPSRDELTKNQVSLTCLVKGFYPSDIAVEWES  
NGQPENNYKTTTPVLDSGDSFFLYSKLTVDKSRWQQGNVFCFSVMHEALHNHYTQKS  
LSLSPGK**

Black: ICOS (a.a. 21-140)

Green: linker

Red: Fc of human IgG1 (a.a.100-330)

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: [\*\*info@bpsbioscience.com\*\*](mailto:info@bpsbioscience.com)

Please visit our website at: [\*\*www.bpsbioscience.com\*\*](http://www.bpsbioscience.com)