

## Data Sheet

### CD40, Fc Fusion (IgG1) Avi-Tag, Biotin Labeled

Human, Recombinant, Fc-Avi-Tag,  
 Biotin Labeled

**Catalog #:** 79102

**Lot #:** 170720

**Conc.:** 1.61 mg/ml

**Formulated in:** 8 mM Phosphate, pH 7.4,  
 110 mM NaCl, 2.2 mM KCl, and 20%  
 glycerol.

**Stability:** At least 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:**  $<1 \text{ EU}/\mu\text{g}$

#### References:

1. Lutgens, E. *et al.*, *Nat. Med.* 1998; **5**:  
 1313-1316.
2. Quezada, S.A., *et al.*, *Immunology.*  
 2003; **22**: 307-328

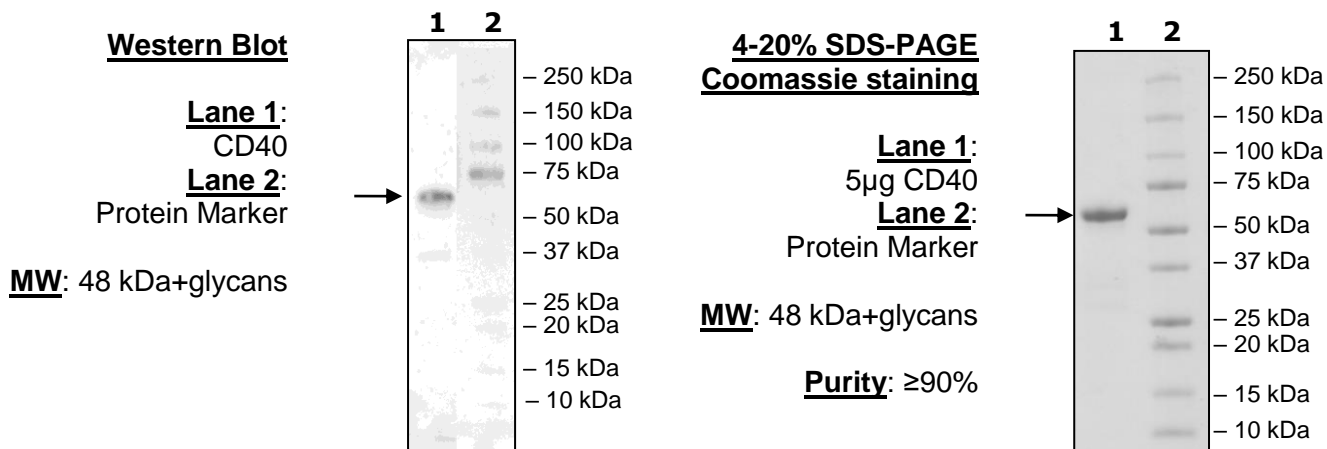
**Description:** Human CD40, also known as  
 Tumor Necrosis Factor Receptor  
 Superfamily Member 5, TNFRSF5, Bp50.  
 GenBank Accession No. NM\_001250, a.a.  
 21-193, fused at the C-terminus with the Fc  
 portion of Human IgG1, with C-terminal Avi-  
 tag™, expressed in a HEK293 cell  
 expression system. MW = 48 kDa. This  
 protein runs at a higher molecular weight  
 due to glycosylation.

**Assay Conditions:** Reaction done  
 according to standard ELISA assay kit  
 protocol (e.g., Catalog # 72005). Buffer is  
 PBS + 0.1% BSA. Blocking buffer is  
 Superblock + 0.05% Tween-20.

Plate was coated with 100 ng/well CD40L  
 (#71191) at  $4^{\circ}\text{C}$  overnight. Various  
 concentrations of biotinylated CD40 (0-1  
 $\mu\text{M}$ ) added for 2h at RT. Strep-HRP added  
 for 1h at RT.

**Applications:** Useful for Western blots,  
 studying protein binding, and for screening  
 small molecules and antibodies.

### Quality Assurance



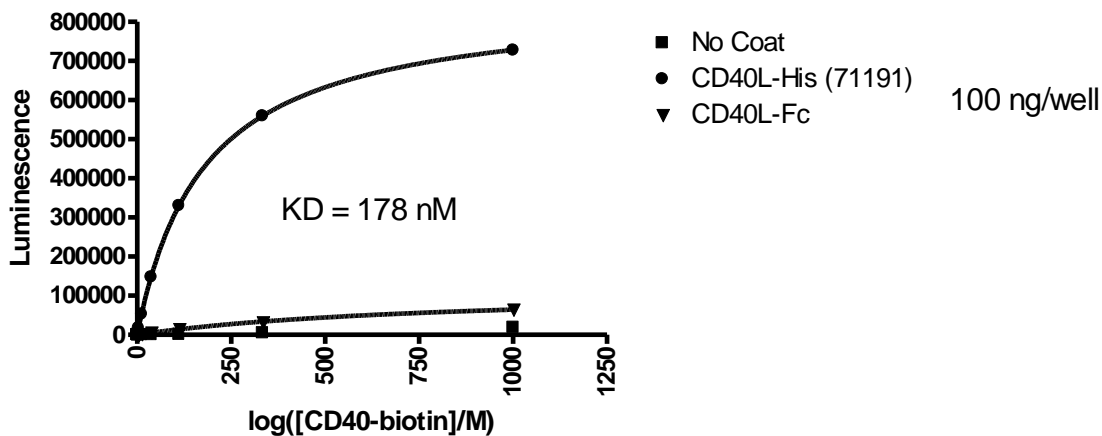
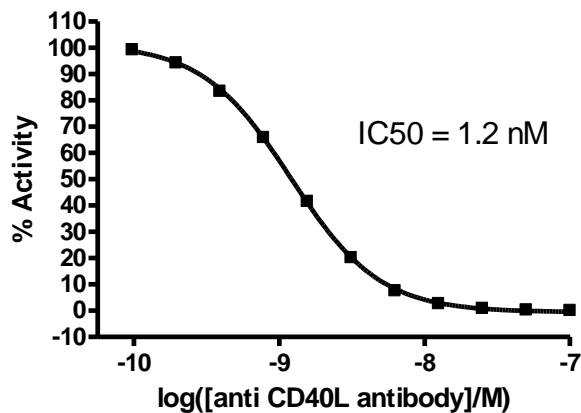
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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)

**CD40: CD40L Interaction**



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)