



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Data Sheet

### CD40L, His-tag

Human, recombinant, N-terminal His-tag

**Catalog #:** 71191

**Lot #:** 191016-1      **Conc.:** 0.87 mg/ml

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

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles.

#### 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 CD40L, also known as CD154, T cell antigen Gp39, TNF related activation protein, TRAP, tumor necrosis factor ligand superfamily member 5, and TNFSF5. GenBank Accession No. NM\_000074, a.a. 116-261 with an N-terminal His-tag, expressed in a HEK293 cell expression system. MW = 18 kDa.

**Assay Conditions:** Assay was performed according to BPS catalog #79257 with CD40 [B] enzyme titrated from 1800 ng/rxn – 0 ng/rxn.

---

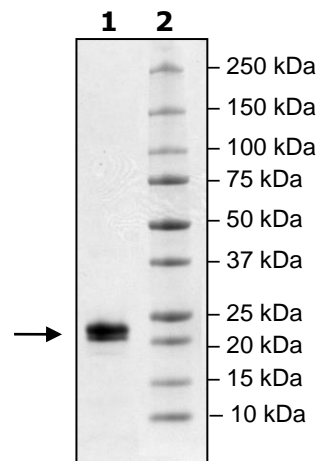
## Quality Assurance

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

**Lane 1:**  
4 $\mu\text{g}$  CD40L  
**Lane 2:**  
Protein Marker

**MW:** 18 kDa

**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 **1.858.202.1401** 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)