

## Data Sheet

### TNFR2, Fc-Fusion (IgG1), His-Avi-Tag Recombinant

Human, Recombinant, Fc fusion protein,  
His-Avi-Tag

**Catalog #:** 79363

**Lot #:** 181106      **Conc.:** 1.29 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.

#### References:

1. Rothe, M., *et al.*, *Cell*. 1995; **83(7)**:  
1243-1252.
2. Xin, C. *et al.*, *Journal of Immunology*.  
2008; **180(10)**: 6467-6471.

#### Description:

Human TNFR2 protein, also known as tumor necrosis factor receptor 2, TNFRSF1B, and CD120b, GenBank Accession No. NM\_001066, a.a. 23-257 with C-terminal His-Avi™-Tag fused to the Fc region of Human IgG1, expressed in a HEK293 cell expression system. MW = 55 kDa.

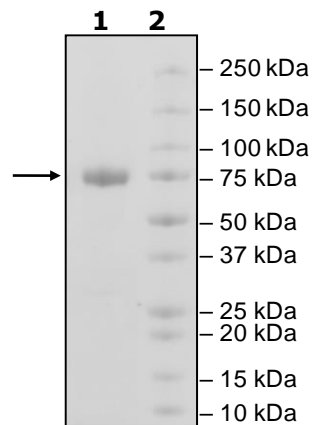
---

### **Quality Assurance**

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

**Lane 1:**  
4 $\mu\text{g}$  TNFR2  
**Lane 2:**  
Protein Marker

**MW:** 55 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: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)