

Data Sheet

TNFR2, Fc-Fusion (IgG1), His-Avi-Tag, Biotin-labeled

Human, Recombinant, Fc fusion protein,
C-terminal His-Avi-Tag, Biotin-labeled

Catalog #: 100205

Lot #: 190321-1 **Conc.:** 2.48 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°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

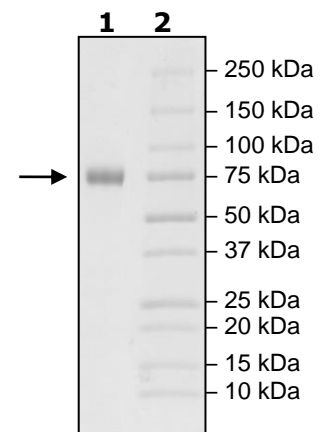
4-20% SDS-PAGE Coomassie staining

Lane 1:
4 μg 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: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com