

Data Sheet

CD30, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled

Human, Recombinant, C-terminal Fc fusion
 Avi-Tag, Biotin-Labeled

Catalog #: 72511

Lot#: 170627-G1 **Conc.:** 0.9 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. Protein may be diluted to
 $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at
 -80°C .

References:

1. Piccinni, M. *et al.*, *Journal of Immunology*. 1995; **1**: 128-133.
2. Smith, C.A. *et al.* *Cell*. 1993; **73(7)**: 1349-1360.

Description:

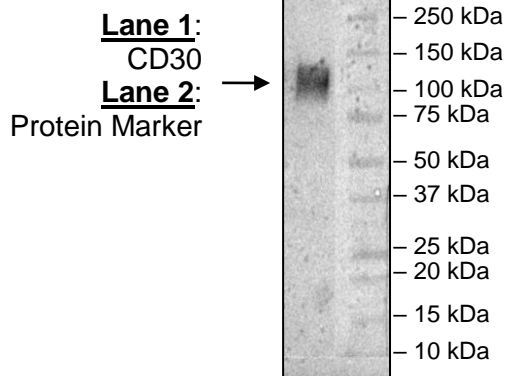
Human CD30, Fc fusion protein, with C-terminal Avi-Tag, Biotin-Labeled, also known as Tumor Necrosis Factor Receptor Superfamily Member 8, and TNFRSF8, GenBank Accession No. NM_001243.4, a.a. 19-379 expressed in a HEK293 cell expression system. MW = 67 kDa. This protein is fused at the C-terminus to the Fc region of human IgG1 and is enzymatically biotinylated using Avi-tag™ technology. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Application:

Useful as a positive control for western blotting.

Quality Assurance

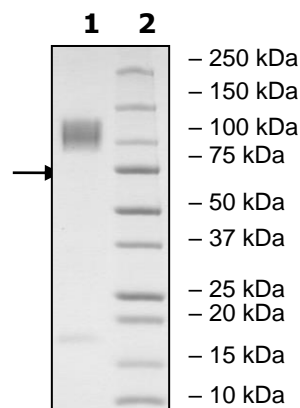
Western Blot



4-20% SDS-PAGE Coomassie staining

Lane 1: CD30-Biotinylated
Lane 2: Protein Marker

MW: 67 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

Please visit our website at: www.bpsbioscience.com