

Data Sheet

CD70 His-Avi-Tag, Biotin-Labeled

Human, Recombinant, C-terminal His-Avi-Tag,
Biotin-Labeled

Catalog #: 100228

Lot#: 181101

Conc.: 0.61 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°C . Avoid
freeze/thaw cycles. Protein may be diluted to
 $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at
 -80°C .

Description:

Human CD70, also known as CD27L,
CD27LG, TNFSF7, TNFSF7G, and CD27
ligand, GenBank Accession No.
NM_001252, a.a. 39-193 (end) with C-
terminal His-Avi-Tag, expressed in a
HEK293 cell expression system. MW = 20
kDa. This protein is enzymatically
biotinylated using Avi-Tag™ technology.
Biotinylation is confirmed to be $>90\%$.

References:

1. Wyzgol, A. *et al.*, *Journal of Immunology*. 2009; **201(10)**: 1851-61.
2. Lorenz, M., *et al.*, *Human Gene Therapy*. 1999; **10(7)**: 1095-103.

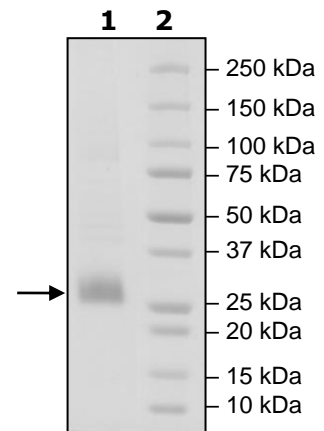
Quality Assurance

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

Lane 1:
4 μg CD27L

Lane 2:
Protein Marker

MW: 20kDa
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