

Data Sheet

CD27(Mouse), Fc-Fusion (Human), Avi-Tag, Biotin

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

Catalog #: 79107

Lot#: 170831-G **Conc.:** 0.91 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. Hendricks, J. *et al.*, *Nat. Immunol.*
2000; **1**: 433-440.
2. Ruprecht, C.R. *et al.*, *J. Exp. Med.*
2005; **201(11)**: 1793-1803.

Description:

Mouse secreted CD27, Fc fusion protein,
also known as Tumor Necrosis Factor
Receptor Superfamily Member 7,
TNFRSF7, and T14, GenBank Accession
No. NM_001033126, a.a. 21-182, with C-
terminal Avi-Tag fused to the Fc portion of
Human IgG1 and expressed in a HEK293
cell expression system. The protein is
enzymatically biotinylated using AviTag™
technology. MW = 47 kDa. This protein runs
at a higher MW by SDS-PAGE due to
glycosylation.

Quality Assurance

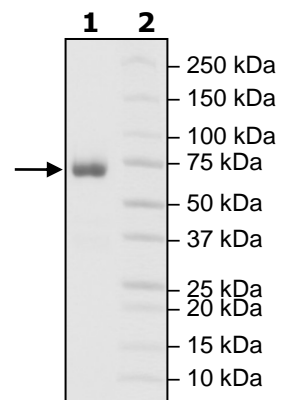
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
Mouse CD27

Lane 2:
Protein Marker

MW: 47 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



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence

TLAPNSCPDKHYWTGGGLCCRMCEPGTFFVKDCEQDRATAAQCDCPCIPGTSFSPDY
HTRPHCESCRHCNSGFLIRNCTVTANAECSCSKNWQCRDQECTECDPPLNPALTR
QPSETPSPQPPPTHLPHGTEKPSWPLHRQLPNSTVYSQRSSHRPLCSSDCIR**DIE**
GRMDPKSCDKTHTCPPCPAPELLGGPSVFLFPPKPKDTLMISRTPEVTCVVDVVS
HEDPEVKFNWYVDGVEVHNAKTKPREEQYNSTYRVVSVLTVLHQDWLNGKEYKCK
VSNKALPAPIEKTISKAKGQPREPQVYTLPPSRDELTKNQVSLTCLVKGFYPSDI
AVEWESNGQPENNYKTTTPVLDSDGSFFLYSKLTVDKSRWQQGNVFCFSVMHEAL
HNHYTQKSLSLSPGKGGGLNDIFEAQKIEWHE

Black = mouse CD27(21-182)
Red = Linker
Blue = Fc of human IgG1 (aa 100-330)
Green = Avi

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