

Data Sheet

CD27, Fc Fusion (IgG1) Avi-Tag, Biotin Labeled

Human, Recombinant, Fc-Avi-Tag,
Biotin Labeled

Catalog #: 100250

Lot #: 181206

Conc.: 1.18 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.

Description: Human CD27, also known as S152, LPFS2, T14, TNFRSF7, and Tp55. GenBank Accession No. NM_001242, a.a. 21-192, fused at the C-terminus of the Fc portion of Human IgG1, with C-terminal Avi-tag™, expressed in a HEK293 cell expression system. MW = 48 kDa. This protein is enzymatically biotinylated using Avi-Tag technology. This protein runs at a higher molecular weight due to glycosylation.

References:

1. Klein, U. *et al.*, *JEM*. 1998; **188(9)**: 1679.
2. Kazunaga, A., *et al.*, *Immunology Today*. 2000; **21(5)**: 204-206

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:

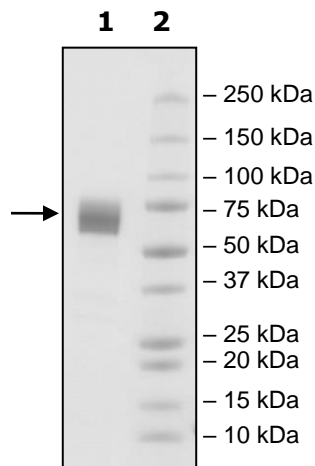
4 μg CD27

Lane 2:

Protein Marker

MW: 48 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 **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com