

Data Sheet

CD27

Human, recombinant, Fc fusion protein

Catalog #: 71176

Lot#: 140528-D **Conc.:** 1.04 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles.

Endotoxin Level: <0.1 EU/ μg

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: Human secreted CD27, Fc fusion protein, also known as Tumor Necrosis Factor Receptor Superfamily Member 7, TNFRSF7, and T14, GenBank Accession No. NM_001242, a.a. 21-192 expressed in a HEK293 cell expression system. MW = 45.8 kDa (monomer). This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Application:

Useful for studying protein binding and screening small molecules.

Quality Assurance

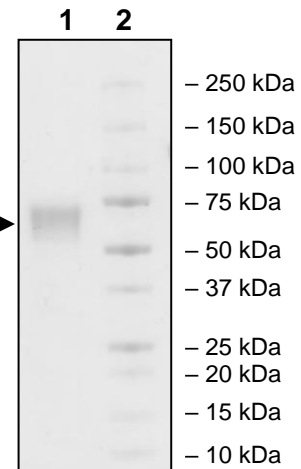
4-20% SDS-PAGE Comassie staining

Lane 1:
3.3 μg CD27

Lane 2: →
Protein Marker

MW: 45.8 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