

Data Sheet

CD277, Fc-Fusion (IgG1) Avi-Tag

Human, Recombinant, Fc fusion protein, with C-terminal Avi-tag™

Catalog #: 100073

Lot #: 180508

Conc.: 0.63 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 ≥ 100 µg/ml in PBS + glycerol and stored at -80°C.

Description:

Human CD227, also known as MUC1, ADMCKD, and MCD, with C-terminal Avi-Tag™ fused to the Fc region of Human IgG1. GenBank Accession No. NM_007048, a.a. 31-254, expressed in a HEK293 cell expression system. MW = 53 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Applications: Useful as a positive control for western blot studies.

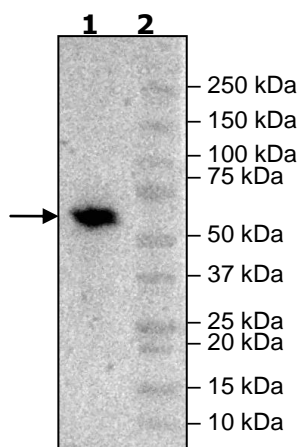
References:

1. Wykes, M., *et al.*, *Journal of Leukocyte Biology*, 2002 Oct; **72(4)**: 692-701.
2. Mukherjee, P., *et al.*, *Journal of Leukocyte Biology*. 2004 Jan; **77(1)**: 90-99.

Quality Assurance

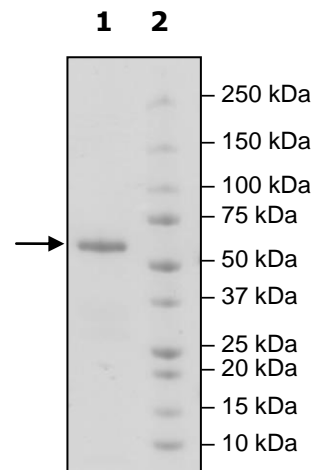
Western Blot

Lane 1:
CD277
Lane 2:
Protein Marker



4-20% SDS-PAGE Coomassie staining

Lane 1:
4µg CD277
Lane 2:
Protein Marker
MW:
53 kDa + glycans
Purity: ≥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