

Data Sheet

CD112, Fc-Fusion, Avi-Tag, Biotin-labeled

Human, Recombinant, C-terminal Fc-Fusion-Avi-Tag, Biotin labeled

Catalog #: 72231

Lot #: 160825

Conc.: 0.77 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. Bottino, C., *et al.*, *J. Exp. Med.* 2003; 198(4): 557-567.
2. Pende, D., *et al.*, *Blood*. 2005; 105(5): 2066-2073.

Description:

Human secreted CD112, also known as Nectin-2, Herpes virus entry mediator B, HVEB, PVRL2, and PRR2, GenBank Accession No. NM_002856, a.a. 32-360 expressed in a HEK293 cell expression system, with C-terminal Fc-Fusion-Avi-tag, this protein is enzymatically biotinylated using Avi-tag™ technology. MW = 64 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

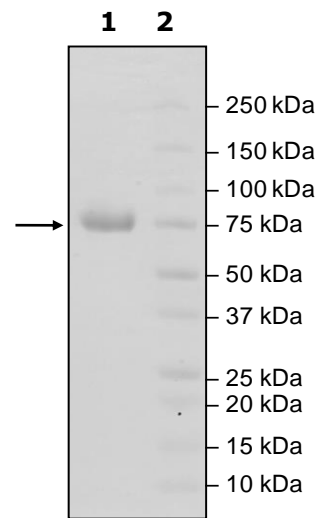
Application:

Useful as a positive control for western blot analysis.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
10 μg CD112
Lane 2:
Protein Marker
MW:
64 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