

Data Sheet

Mouse CD137, Fc fusion (mlgG2a), Biotin-labeled

Mouse, recombinant, Fc fusion protein with C-terminal AviTag™

Catalog #: 71255

Lot #: 160309

Conc.: 1.27 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.

References:

1. Wilcox, R.A., *et al.*, *J. Clin. Invest.* 2002; **109(5)**: 651-659.
2. Foell, J. *et al.*, *J. Clin. Invest.* 2003; **111(10)**: 1505-1518.

Description:

Mouse secreted CD137, also known as Tumor Necrosis Factor Receptor Superfamily member 9, TNFRSF9, and ILA, GenBank Accession No. NM_001561, a.a. 24-187, fused at the C-terminus to the Fc portion of mouse IgG2a, with a C-terminal AviTag™ expressed in a HEK293 cell expression system. This protein is enzymatically biotinylated using AviTag™ technology. MW = 46.7 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.

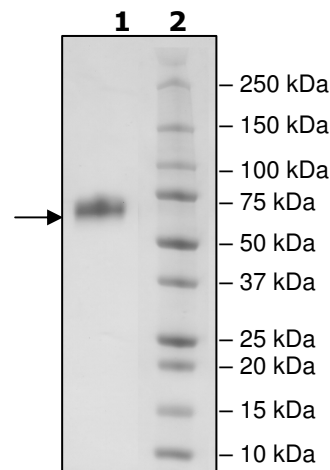
Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
3 µg mCD137

Lane 2:
Protein Marker

MW: 46.7 kDa
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**

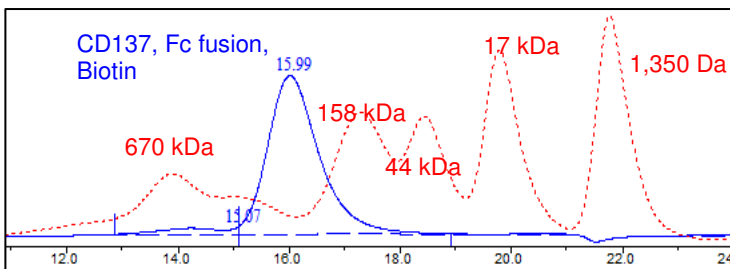
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Gel Filtration Curve



<10% aggregation

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**](mailto:info@bpsbioscience.com)
Please visit our website at: [**www.bpsbioscience.com**](http://www.bpsbioscience.com)