

Data Sheet

CD137, Fc Fusion, Avi-tag **Biotin-Labeled**

Human, Recombinant protein,
 C-terminal Fc-Avi-tag

Catalog #: 71171

Lot #: 191106-G **Conc.:** 0.91 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.

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.

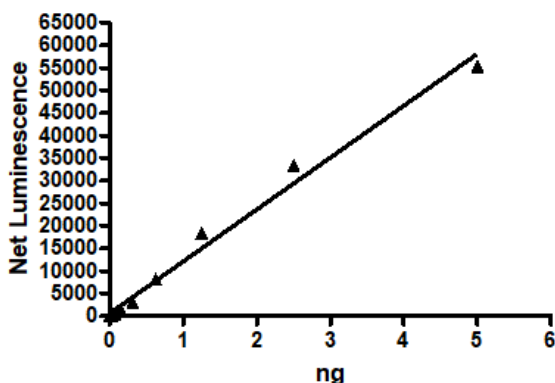
Description: Human secreted CD137, also known as Tumor Necrosis Factor Receptor Superfamily member 9, TNFRSF9, and ILA, GenBank Accession No. NM_001561, a.a. 24-186, fused at the C-terminus to the Fc portion of human IgG1, expressed in a HEK293 cell expression system and enzymatically biotin-labeled using Avi-tag™ technology. MW = 46 kDa (monomer). This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Assay Conditions: Assay was performed according to BPS catalog #72025 with CD137[B] enzyme titrated from 40 ng/rxn – 0 ng/rxn.

Application: Useful for studying protein binding and screening small molecules.

Quality Assurance

CD137L:CD137[B] Interaction

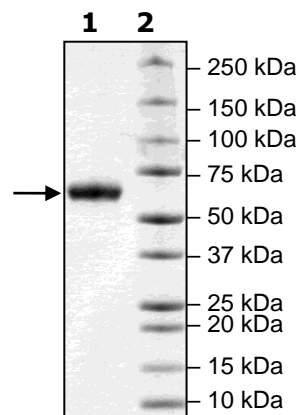


4-20% SDS-PAGE **Coomassie staining**

Lane 1:
 2 µg Cd137
Lane 2:
 Protein Marker

MW: 46 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 **1.858.202.1401** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com