

## Data Sheet

### Mouse CD137, Fc fusion (hlgG1)

Mouse, recombinant, Fc fusion protein

**Catalog #:** 71250

**Lot #:** 160217

**Conc.:** 2.0 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  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Storing diluted protein is not recommended, if necessary, include carrier protein (BSA 0.1 – 0.5%).

**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\_011612, a.a. 24-187 fused at the C-terminus to the Fc portion of human IgG1, expressed in a HEK293 cell expression system. MW = 44 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.

---

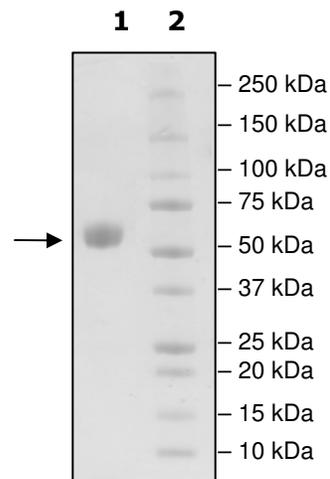
### Quality Assurance

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

**Lane 1:**  
4  $\mu\text{g}$  mCD137

**Lane 2:**  
Protein Marker

**MW:** 44 kDa  
**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**

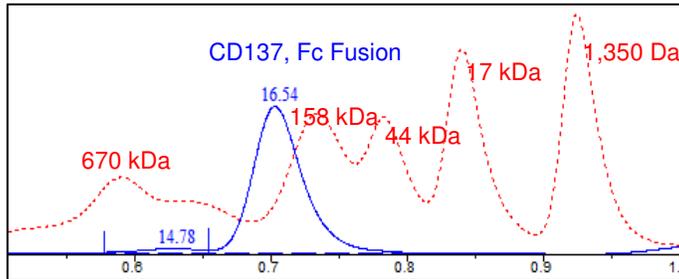
Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://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 Data



<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)