

Data Sheet

OX40 (CD134), Fc-Fusion

Human, Recombinant, Fc fusion protein

Catalog #: 71175

Lot #: 190612 **Conc.:** 1.30 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. Rogers, P.R. *et al.*, *Immunity*. 2001; **15(3)**: 445-455.
2. Ito, T. *et al.*, *J. Exp. Med.* 2005; **202(9)**: 1213-1223.

Description:

Human secreted OX40, Fc fusion protein, also known as Tumor Necrosis Factor Receptor Superfamily Member 4, TNFRSF4, and CD134, GenBank Accession No. NM_003327, a.a. 29-216 expressed in a HEK293 cell expression system. MW = 47 kDa (monomer). This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Quality Assurance

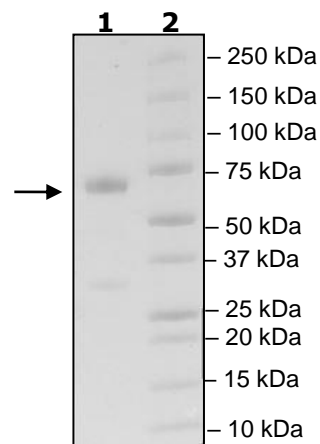
4-20% SDS-PAGE Coomassie staining

Lane 1:
OX40

Lane 2:
Protein Marker

MW: 47 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 **1.858.202.1401** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com