

Data Sheet

OX40L, His-Avi-Tag

Human, Recombinant, N-terminal His-Avi-tag

Catalog #: 100808

Lot #: 200715

Conc.: 1.07 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.

Description: Human OX40L, also known as tumor necrosis factor ligand superfamily member 4 (TNFSF4) and CD252, GenBank Accession No. NM_003326, a.a. 51-183(end) with a N-terminal His-Avi-tag, expressed in a HEK293 cell expression system. MW = 23 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

References:

1. Linton, P.J., *et al.*, *J. Exp. Med.* 2003; **197(7)**: 875-883.
2. Arestides, R.S.S., *et al.*, *Eur. J. Immunol.* 2002; **32(10)**: 2874-2880.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:

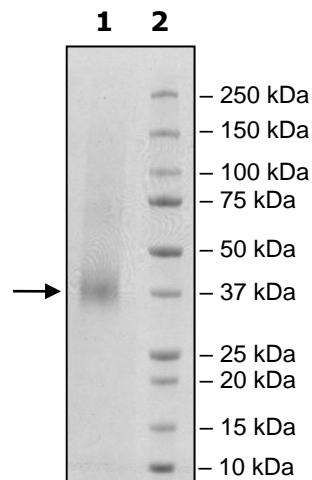
4 μg OX40L

Lane 2:

Protein Marker

MW: 23 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