

Data Sheet

Human LTBR, Fc Fusion HiP™ Recombinant

Human, recombinant, Fc fusion protein

Catalog #: 71112

Lot#: 131113-D Conc.: 0.70 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, and 20%
glycerol

Stability: >6 months at -80°C. *Avoid
freeze/thaw cycles.*

References:

1. Crowe, P.D., *et al.*, *Science*. 1994
Apr 29;**264(5159)**:707-10.
2. Wu, M.Y., *et al.*, *J Biol Chem*. 1999
Apr 23;**274(17)**:11868-73.

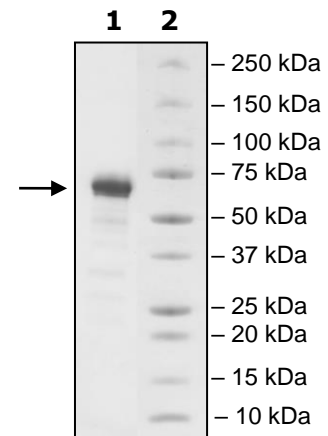
Description:

Human secreted lymphotoxin beta receptor (LTBR)-Fc fusion protein, also known as, Tumor Necrosis Factor C Receptor (TNFCR), and CD18, GenBank Accession No. NM_002342, a.a 28-219, expressed in a HEK293 cell expression system. MW = 48 kDa (monomer). This protein runs at a higher apparent M.W. by SDS-PAGE due to glycosylation.

Application:

Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

**4-20% SDS-PAGE
Coomassie staining****Lane 1:**
3.5 µg LTBR-Fc**Lane 2:**
Protein Marker**MW:** 48 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 **1.858.202.1401** Fax **1.858.481.8694**Or you can Email us at: support@bpsbioscience.comPlease visit our website at: www.bpsbioscience.com