

## Data Sheet

### Mouse LTBR, Fc Fusion HiP™ Recombinant

Mouse, Recombinant, Fc fusion protein

**Catalog #:** 71122**Lot #:** 140318-2      **Conc.:** 2.76 mg/ml**Formulated in:** 10 mM phosphate, 138 mM NaCl, 2.7 mM KCl, pH 7.4.**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . *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.

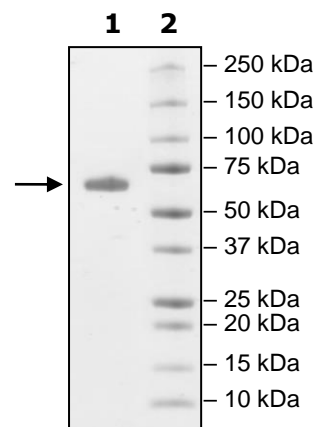
**Description:**

Mouse secreted lymphotoxin beta receptor (mLTBR)-Fc fusion protein, also known as, Tumor Necrosis Factor C Receptor (TNFCR), and CD18, GenBank Accession No. NM\_010736, a.a 28-221, expressed in a HEK293 cell expression system. MW = 48.8 kDa (monomer). Endotoxin level  $\sim 0.0007$  EU/ $\mu\text{g}$ . This protein runs at a higher apparent M.W. by SDS-PAGE due to glycosylation.

**Application:**

Useful for the study of protein-protein interaction and in vitro and in vivo protein function studies.

## Quality Assurance

**4-20% SDS-PAGE  
Coomassie staining****Lane 1:**  
2  $\mu\text{g}$  TBR-Fc**Lane 2:**  
Protein Marker**MW:** 48.8 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 **1.858.202.1401** Fax **1.858.481.8694**Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)