

Data Sheet

BTNL8, Fc-fusion (IgG1), Avi-Tag

Human, recombinant, C-terminal Fc-Avi-Tag,
Catalog #: 100467

Lot #: 190828 **Conc.:** 0.59 mg/ml

Formulation: 8 mM Phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol.

Stability: Stable for at least 6 months at -80°C. Avoid freeze/thaw cycles.

References:

1. Chapoval, A.I., *et al.*, *Molecular Immunology*. 2013 Dec; **56(4)**: 819-828.
2. Riquelme, W.L.. *et al.*, *Transplantation*. 2014 Jul; **98**: 396.

Description: Human BTNL8, also known as Butyrophilin-like protein 8, GenBank Accession No. NM_024850, a.a. 18-238, fused to the C-terminus at the Fc portion of human IgG1 followed by a C-terminal Avi-tag™ and expressed in a HEK293 cell expression system. MW=54 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Quality Assurance

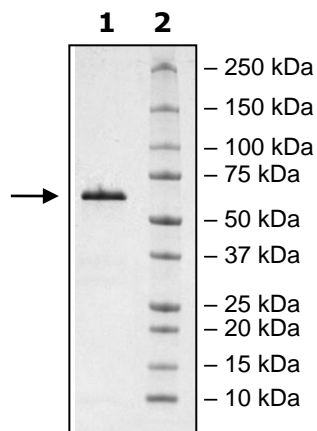
4-20% SDS-PAGE
Coomassie staining

Lane 1:
 4µg BTNL8

Lane 2:
 Protein Marker

MW: 54 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