

Data Sheet

BTN1A1 Avi-His-Tag

Human, Recombinant, C-terminal Avi-His-Tag

Catalog #: 100469

Lot#: 191125

Conc.: 1.0 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 BTN1A1, also known as Butyrophilin subfamily 1 member A1, GenBank Accession No. NM_001732, a.a. 27-242 with a C-terminal Avi-His-Tag™, expressed in a HEK293 cell expression system. MW=27 kDa. This protein runs at a higher molecular weight due to glycosylation.

References:

1. Ogg, S.L. *et al.*, *PNAS*. 2004; **101(27)**: 10084-10089.
2. Smith, I. A. *et al.*, *J. Immunology*. 2010; **184(7)**: 3514-3525.

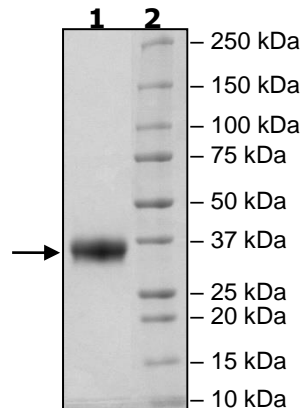
Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg BTN1A1
Lane 2:
Protein Marker

MW: 27 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: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com