

Data Sheet

TIM-3, Avi-His-Tag

Human, Recombinant, C-terminal Avi-His-Tag
Catalog#: 71262
Lot#: 150729-2 **Conc.:** 0.55 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. Protein may be diluted to $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at -80°C .*

References:

1. Zhu, C. *et al.*, *Nat. Immunol.* 2005;
6:1245-1252.
2. Anderson, A.C. *et al.*, *Science.* 2007;
318(5853):1141-1143.

Description:

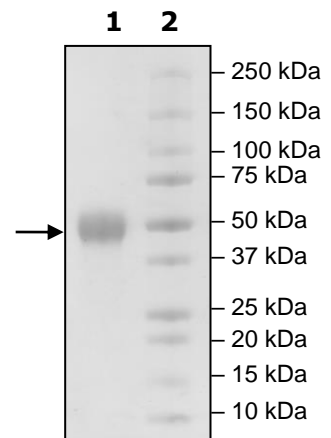
Human TIM-3, also known as T-cell immunoglobulin mucin receptor 3, T-cell membrane protein 3, TIMD-3, Hepatitis A virus cellular receptor 2, HAVCR-2, and KIM-3. GenBank Accession No. NM_032782, a.a. 22-200 with a C-terminal Avi-His-tag, expressed in a HEK293 cell expression system. MW = 23 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Quality Assurance

**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
3 μg Tim3
Lane 2:
Protein Marker

MW: 23 kDa+glycans
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 **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com