

Data Sheet

TIM-3, Biotin-labeled, His-tag

Human, recombinant, C-terminal Avi-biotin, His- tag

Catalog#: 71237

Lot#: 161025

Conc.: 0.62 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:

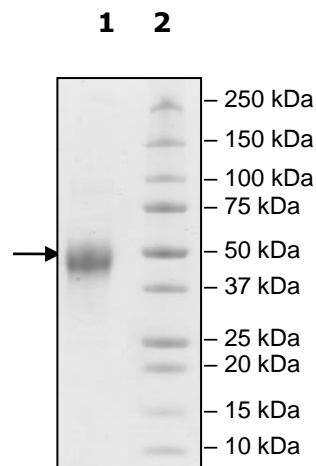
Human TIM-3, also known as T-cell immunoglobulin mucin receptor 3, T-cell membrane protein 3, TIMD-3, Hepatitis A virus cellular receptor 2, HAVCR-2, and KIM-3. GenBank Accession No. NM_032782, a.a. 22-200 with a C-terminal Avi-biotin, His-tag, expressed in a HEK293 cell expression system. MW = 23.3 kDa. This protein is enzymatically biotinylated using Avi-tag™ technology. Biotinylation is confirmed to be greater than 90%. 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: 23.3kDa
Purity: $\geq 90\%$



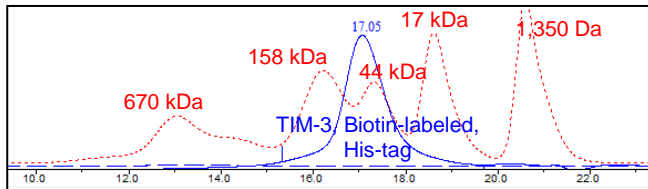
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Gel Filtration Curve



<5% aggregation

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