

Data Sheet

TIM-3, Avi-Tag, Biotin Labeled

Human, Recombinant, C-terminal Avi-Tag
Catalog#: 79096
Lot#: 170811 **Conc.:** 0.23 mg/ml

Formulated in: 40mM Tris pH8.0, 110mM NaCl, 2.2mM 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:

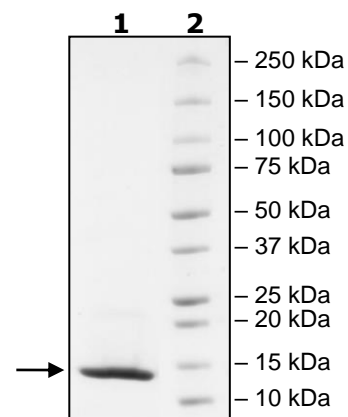
Human TIM-3, also known as T-cell immunoglobulin mucin receptor 3, T-cell membrane protein 3, TIMD-3, Hepatitis A virus cellular receptor 2, HAVCR-2, and KIM-3. GenBank Accession No. NM_032782, a.a. 24-130 with a C-terminal Avi-tag, expressed in an E. coli cell expression system and enzymatically biotin-labeled using AviTag™ technology. MW = 14 kDa.

Quality Assurance

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

Lane 1:
6 μg TIM3
Lane 2:
Protein Marker

MW: 14 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 **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com