

Data Sheet

TIM-3, FLAG-Tag

Human, Recombinant, C-terminal FLAG- tag
Catalog#: 71180
Lot#: 170531 **Conc.:** 0.07 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 .*

Endotoxin Level: $<1 \text{ EU}/\mu\text{g}$

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 with a C-terminal FLAG-tag. GenBank Accession No. NM_032782.4, a.a. 22-200 expressed in a HEK293 cell expression system. MW = 21 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:

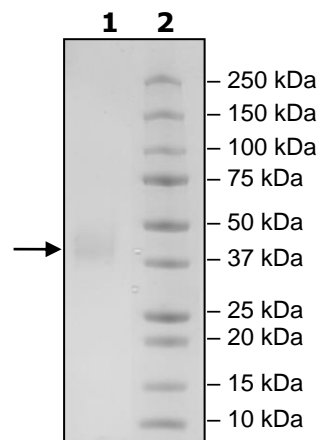
2 μg Tim3

Lane 2:

Protein Marker

MW: 21 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



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence

SEVEYRAEVLGQNAAYLPCFYTPAAPGNLVPVCWGKGACPVFECGNVVLRTDERDVN
YWTSRYWLNGLDFRKGDVSLTIENVTLADSGIYCCRIQIPGIMNDEKFNLKLVIKP
AKVTPAPTRQRDFTAAFPRLTTRGHGPAETQTLGSLPDINLTQISTLANELRDS
RLANDLRDSGATIRDYKDDDDK

Black: TIM-3 (a.a. 22-200)

Blue: FLAG

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**](mailto:info@bpsbioscience.com)
Please visit our website at: [**www.bpsbioscience.com**](http://www.bpsbioscience.com)