

Data Sheet

B7-H3, Avi-His-Tag, Biotin-Labeled

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

Catalog #: 79367

Lot#: 171109 **Conc.:** 1.9 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:

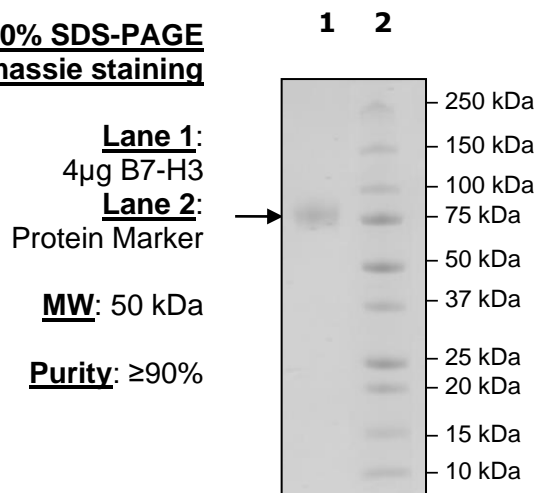
1. Chapoval, A., et al., Nature Immunology. 2001; 2(3): 269-74.
2. Sun, M., et al., J. Immunol. 2002; 168(12):6294-7.

Description:

Human secreted B7-H3, known as Cluster of Differentiation 276 and CD276, GenBank Accession No. NM_001024736, a.a. 29-466, with C-terminal Avi-His-Tag expressed in a HEK293 expression system. This protein is enzymatically biotinylated using AviTag™ technology. MW = 50 kDa.

Quality Assurance

4-20% SDS-PAGE Coomassie staining



BPS BIOSCIENCE 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**

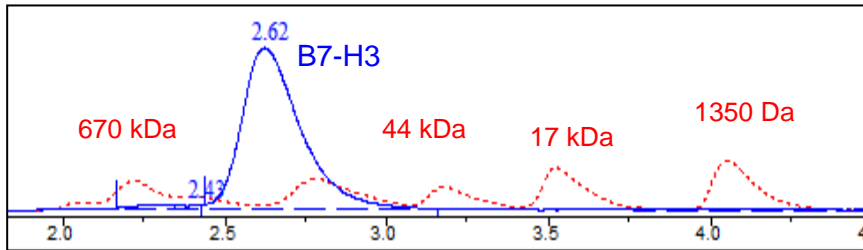
Or you can Email us at: info@bpsbioscience.com

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

Gel Filtration Curve



<10% Aggregation

BPS BIOSCIENCE 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