

Data Sheet

B7-H4, His-tag

Human, recombinant, C-terminal His-tag
Catalog #: 71144
Lot#: 140331 Conc.: 1.1 mg/ml

Formulated in: 40 mM Tris, pH8.0, 110 mM NaCl, 2.4 mM KCl, and 20% glycerol.

Stability: >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: <0.1 EU/ μg

References:

1. Byrum, J.N., *et al.*, *J. Immunol.* 2013 Sep 15;**191(6)**:3073-3081.
2. Legat, A., *et al.*, *Front Immunol.* 2013 Dec 19;**4**:455.

Description:

Human secreted B7-H4, also known as B7x and B7S1, with C-terminal His-tag GenBank Accession No. NM_024626, a.a. 29-258 expressed in a HEK293 cell expression system MW = 26.7 kDa (monomer). This protein runs at a higher M.W. by SDS-PAGE due to glycosylation.

Application:

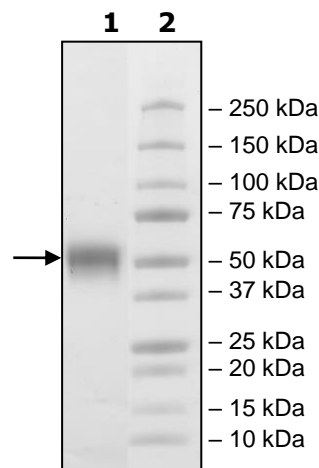
Useful for the study of protein-protein interaction and in vitro and in vivo protein function studies.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
2.8 μg B7-H4
Lane 2:
Protein Marker

MW: 26.7 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



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

Sequence

**FGISGRHSITVTTVASAGNIGEDGILSCTFEPDIKLSDIVIQWLKEGVLGLVHEFKEGKD
ELSEQDEMFRGRTAVFADQVIVGNASRLKKNVQLTDAGTYKCYIITSKGKGNANLEYK
TGAFSMPEVNVVDYNASSETLRCEAPRWFPQPTVVWASQVDQGANSFSEVSNTSFELN
SENVTMKVVSVLNVNTINNTYSCMIENDIAKATGDIKVTSEIKRRSHLQLLNSKAHHH
HHHHHHH**

Black: B7H4 (a.a. 29-258)

Red: His (10x)

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)