

Data Sheet

B7-1 (CD80), Fc fusion, Avi-Tag, Biotin-labeled

Human, Recombinant, Fc fusion protein,
C-terminal Avi-Tag, Biotin labeled.

Catalog #: 71114-2

Lot #: 160901 **Conc.:** 0.86 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:

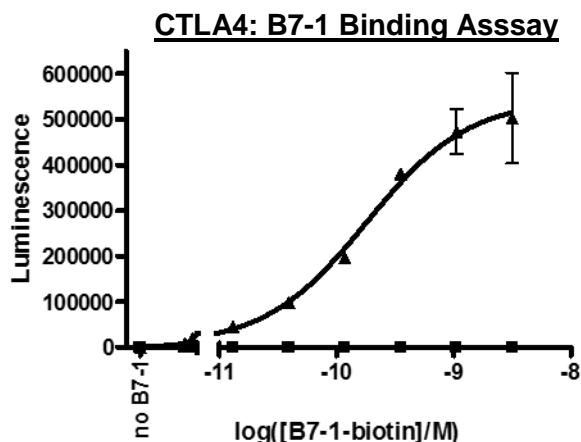
- Byrum, J.N., Van Komen, J.S.,
Rodgers, W., *J. Immunol.* 2013 Sep
15;191(6):3073-3081.
- Legat, A., et al., *Front Immunol.* 2013
Dec 19;4:455.

Description: Human secreted B7-1-Fc
fusion protein, also known as CD80, with C-
terminal Avi-Tag. GenBank Accession No.
NM_005191, a.a 35-242 expressed in a
HEK293 cell expression system and
enzymatically biotin-labeled using Avi-tag™
technology. MW = 53 kDa (monomer). This
protein runs at a higher M.W. by SDS-
PAGE due to glycosylation.

Assay Conditions: Reaction done
according to CTLA4:B7-1[biotin] assay kit
(Catalog # 72009) protocol. Buffer is PBS +
0.1% BSA. Blocking buffer is Superblock +
0.05% Tween-20. Coat protein was added
ON at 4 deg 100 ng/well. Various
concentrations of biotinylated ligand (0-13.5
ng/ul, 0.5 ng/ul is the recommended kit
condition) added for 2h at RT. Strep-HRP
added for 1h at RT.

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

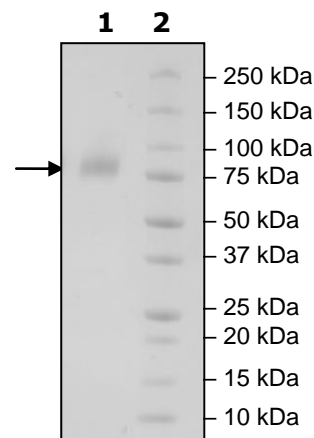
Quality Assurance



4-20% SDS-PAGE Coomassie staining

Lane 1:
3 μg B7-1
Lane 2:
Protein Marker

MW: 53 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**

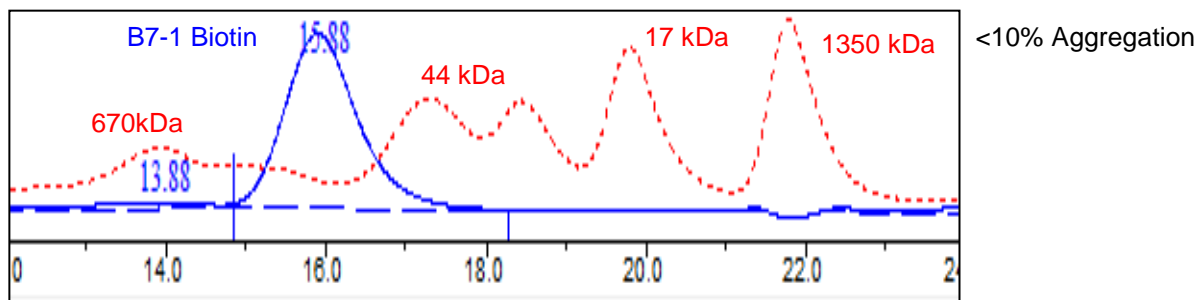
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



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



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