

Data Sheet

B7-1 (CD80), Fc Fusion

Human, recombinant, Fc fusion protein

Catalog #: 71125

Lot #: 150310

Conc.: 2.29 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 $\geq 100\text{ }\mu\text{g/ml}$ in PBS + glycerol and stored at -80°C .

References:

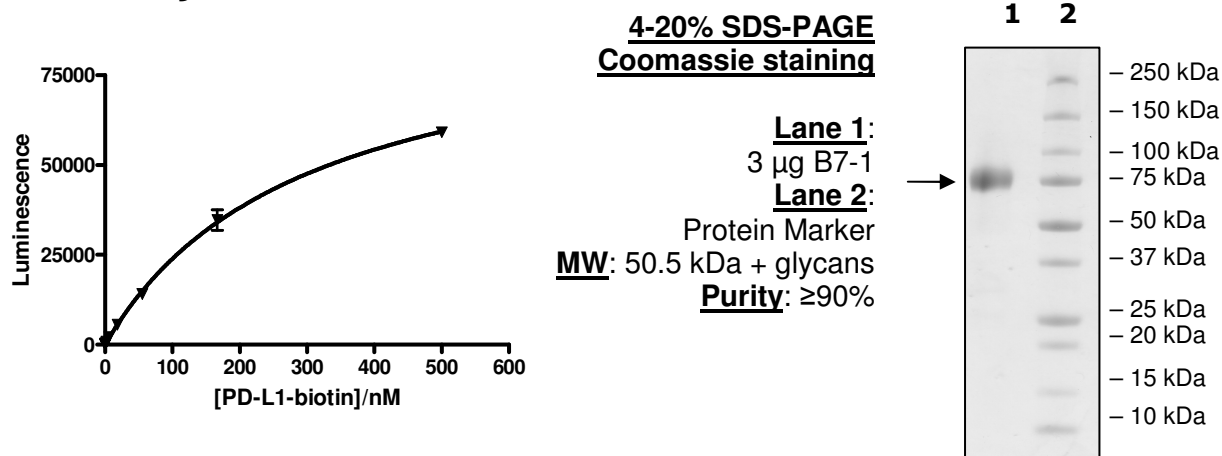
1. Byrum, J.N., *et al.*, *J. Immunol.* 2013; **191(6)**: 3073-3081
2. Legat, A., *et al.*, *Front. Immunol.* 2013; **19(4)**: 455.

Description: Human secreted B7-1-Fc fusion protein, also known as CD80. GenBank Accession No. NM_005191, a.a. 35-242 expressed in a HEK293 cell expression system. MW = 50.5 kDa (monomer). This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Assay Description: B7-1 binding to PD-L1 was determined by coating 100 ng of B7-1 per well and adding the indicated amount of PD-L1-biotin (BPS Cat. #71105) and incubating for 2 hr at RT. Strep-HRP was added and allowed to incubate at room 1 hr at RT, followed by HPR chemiluminescent substrate addition and plate reading.

Application: Useful for the study of protein binding studies and screening inhibitors.

Quality Assurance



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**

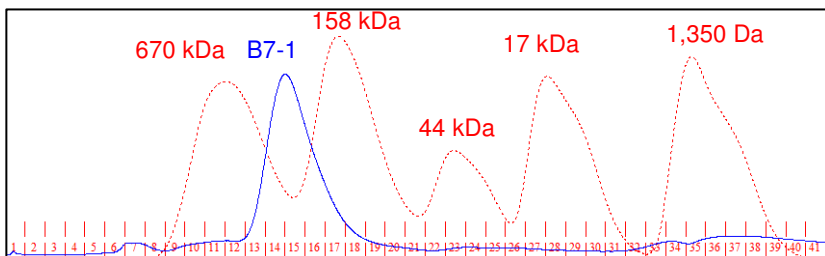
Or you can Email us at: info@bpsbioscience.com

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

Gel Filtration Curve:



<5% Aggregation

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)