



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

Data Sheet

Anti-PD-L1 Antibody, Biotin-labeled

Human, Recombinant, Monoclonal Anti-PD-L1
Catalog #: 71214
Lot #: 141027-BDS **Conc.:** 1.07 mg/ml

Host Species | Isotype:

Human | IgG1

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, and 20%
glycerol.

Endotoxin Level: <1 EU/μg

Stability: Stable for at least 12 months at
-80°C. *Avoid freeze/thaw cycles. Store
diluted antibody with carrier protein (BSA 0.1
– 0.5%) at 4°C, stable for 2 weeks.*

References:

1. Freeman, G.J., *et al.*, *J. Exp. Med.* 2000;
192(7): 1027-1034
2. Dong, H., *et al.*, *Nat. Med.* 1999; **5(12):**
1365-1369.

Description: Purified, recombinant,
humanized anti-PD-L1 (IgG1) antibody with
C-terminal Avi-tag, expressed in HEK293
cells. Purified antibody was enzymatically
biotinylated using Avitag™ technology. This
antibody reacts with human PD-L1 and has
not been tested for cross-reactivity with other
species.

Purification: Protein A affinity
chromatography from HEK293 supernatants.

Applications: Western blot, ELISA

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