

Data Sheet

Anti-PD-1

Recombinant, Murine/Human Monoclonal
Anti-PD-1

Catalog #: 71120

Lot #: 160420 **Conc.:** 1.47 mg/ml

Host Species | Isotype:

Murine/Human | IgG1k

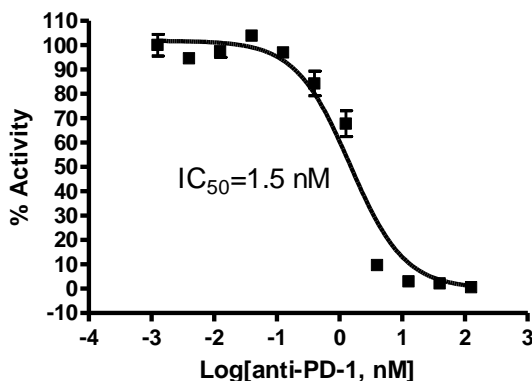
Formulation: Sterile filtered in 8 mM phosphate pH7.4 110 mM NaCl, 2.2 mM KCl, and 20% glycerol.

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. Shinohara, T., *et al.*, *Genomics*. 1994 Oct;**23(3)**:704-6.
2. James, E.S., *et al.*, *Genes Immun*. 2005 Aug;**6(5)**:430-7.

Image(s):



Description: Neutralizing recombinant murine/human chimeric antibody recognizing the PD-L1/PD-L2 binding region of human PD-1. This antibody does not cross-react with mouse PD-1, but does cross-react with monkey (*M. fascicularis*) PD-1. It has not been tested with other species.

IC50: 1.5 nM using the PD-1[Biotinylated]:PD-L1 Inhibitor Screening Assay Kit (BPS Cat. #72005).

Assay Conditions: Antibody was added to PD-1[Biotinylated]:PD-L1 or PD-1[Biotinylated]:PD-L2 binding assay at various concentrations. Assay was performed according to the protocol in BPS Cat. #72005 or #72006, respectively.

Purification: Protein A affinity chromatography from HEK293 supernatants.

Application: Useful for inhibition of PD-1:PD-L1 and PD-1:PD-L2 interactions

Figure 1. Wells were coated with 100 ng of PD-L1 (BPS Cat. #71104). 10 ng of PD-1, Biotin-labeled (BPS Cat. #71109) and varying concentrations of PD-1 Neutralizing Antibody (BPS Cat. #71120) were added and plate was incubated at room temperature for 2 hours. Straptavidin-HRP was added, followed by HRP substrate and chemiluminescence was read.

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

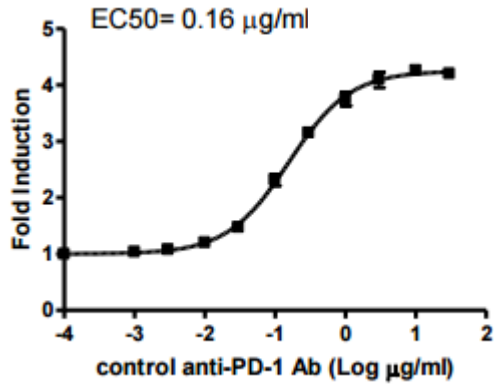


Figure 2. PD-1 antibody was confirmed to neutralize PD-1 mediated NFAT suppression using BPS PD-1 Reporter Cell Line (BPS Cat. #60535)

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