

Data Sheet

Anti-PD-1 Neutralizing Antibody, PE-labeled

Recombinant, Human/Murine Monoclonal
 Anti-PD-1

Catalog #: 71290-2

Lot #: 160323

Conc.: 5.14 mg/ml

Host Species | Isotype:

Murine/Human | human IgG1k

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

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

References:

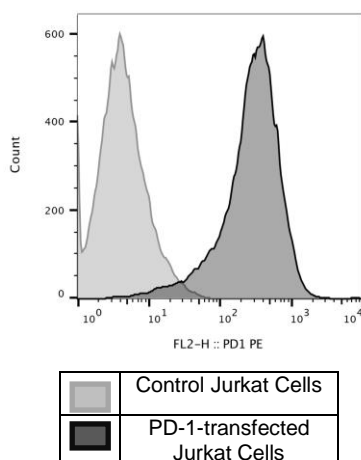
1. Shinohara, T., *et al.*, *Genomics*. 1994
 Oct;**23**(3):704-6.
2. James, E.S., *et al.*, *Genes Immun*. 2005
 Aug;**6**(5):430-7.

Description: Phycoerythrin-labeled anti-PD-1
 neutralizing recombinant human/murine
 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.

Recommended Assay Conditions: For flow
 cytometry, we recommend incubating cells on
 ice for 30 minutes with 1-20 µg/ml of anti-
 human PD-1 antibody, PE-labeled.

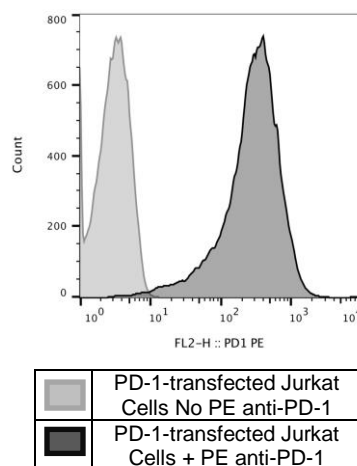
Purification: Protein A affinity chromato-
 graphy from HEK293 supernatants.

Application: Useful for labeling cells
 expressing PD-1 for flow cytometry and
 immunofluorescence microscopy.



Control Jurkat cells and PD-1-
 expressing stable Jurkat cells
 (BPS Cat. #60535) were
 treated with PE anti-human
 PD-1 antibody (left).

PD-1-expressing stable Jurkat
 cells (BPS Cat. #60535) with
 or without PE Anti-human PD-
 1 antibody (right).



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