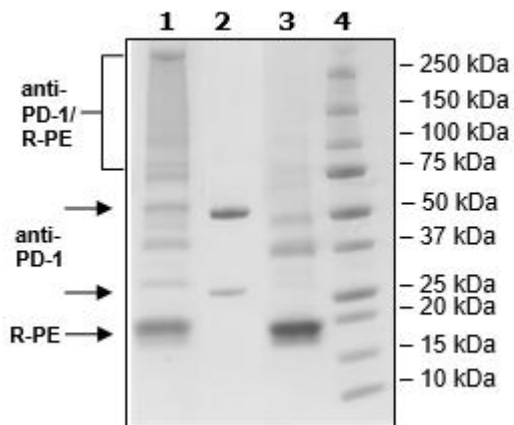


## Product Information

<b>Description:</b>	Phycoerythrin-labeled anti-PD-1 neutralizing recombinant human/murine chimeric antibody recognizing the PD-L1/PD-L2 binding region of human PD-1.
<b>Label:</b>	R-Phycoerythrin (PE) is an oligomeric protein complex (270 kDa) from red algae that exhibits intensely bright red-orange fluorescence with high quantum yields. The complex consists of six heterodimers, $\alpha$ subunit (18 kDa) and $\beta$ -subunit (20 kDa), and an additional $\gamma$ -subunit (34 kDa). PE is covalently attached randomly through lysines on the target protein.
<b>Concentration:</b>	5.14 mg/ml
<b>Host Species:</b>	Murine/Human
<b>Isotype:</b>	IgG1 $\kappa$
<b>Clonality:</b>	Monoclonal
<b>Cross Reactivity:</b>	This antibody does not cross-react with mouse PD-1, but does cross-react with monkey ( <i>M. fascicularis</i> ) PD-1. It has not been tested with other species.
<b>Formulated In:</b>	8 mM phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
<b>Expression System:</b>	HEK293
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 12 months at -80°C. Avoid freeze/thaw cycles. Protect from light.
<b>Storage:</b>	-80°C
<b>MW:</b>	Heavy Chain: 49 kDa + glycans + PE, Light Chain: 23 kDa + PE
<b>Purity:</b>	≥90%
<b>Purification:</b>	Protein A affinity chromatography from HEK293 supernatants.
<b>Assay Conditions:</b>	For flow cytometry, we recommend incubating cells on ice for 30 minutes with 1-20 $\mu$ g/ml of anti-human PD-1 antibody, PE-labeled.
<b>Applications:</b>	Useful for labeling cells expressing PD-1 for flow cytometry and immunofluorescence microscopy.

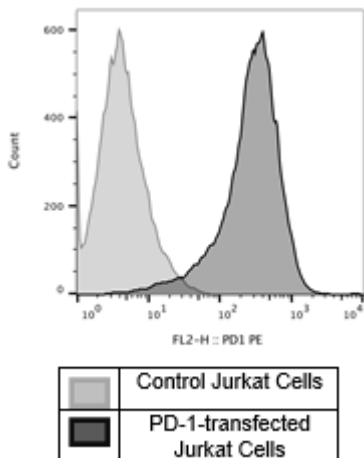
## Quality Control Data

### 4-20% SDS-Page Coomassie Staining

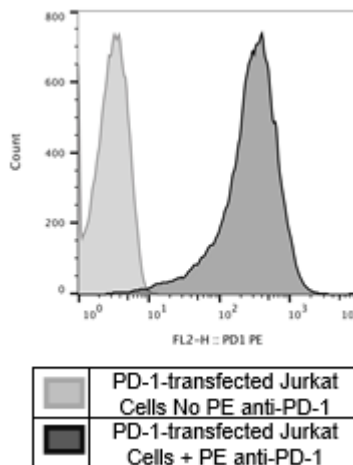


**Lane 1:** 8 µg anti-PD-1/ R-PE  
**Lane 2:** 4 µg anti-PD-1  
**Lane 3:** 4 µg R-PE  
**Lane 4:** Protein Marker

### FACS Assay



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



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