

## Data Sheet

### Anti-GITR, PE-labeled

Recombinant, Human Monoclonal  
 Anti-GITR

**Catalog #:** 71295-2

**Lot #:** 160812

**Conc.:** 1.68 mg/ml

**Host Species | Isotype:**

Human | Human IgG1

**Formulated in:** 8 mM Na Phosphate pH 7.4,  
 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:**

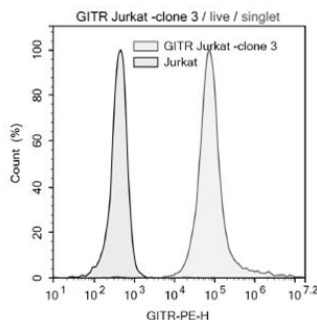
1. Cohen, A., *et al.*, *Cancer Research*.  
 2006 May; **66(9)**: 4904-4912.
2. Scumpia, P.O., *et al.*, *Blood*. 2007 Oct;  
**110(10)**:3673-3681.

**Description:** Phycoerythrin-labeled anti-GITR  
 recombinant human antibody recognizing the  
 GITRL binding region of human GITR. This  
 antibody has not been tested for cross-reactivity  
 with any other species.

**Recommended Assay Conditions:** For flow  
 cytometry, we recommend incubating cells  
 with 20 µg/ml of anti-human PD-1 antibody,  
 PE-labeled for one hour at 4°C, followed by  
 two washes with FACS buffer.

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

**Application:** Useful for labeling cells  
 expressing GITR for flow cytometry and  
 immunofluorescence microscopy.



	GITR Jurkat – clone 3
	Control Jurkat Cells

Control Jurkat cells and GITR expressing stable Jurkat cells were treated with PE anti-human GITR antibody and analyzed for fluorescence by flow cytometry.

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)