

Data Sheet

Anti-GITR, PE-labeled

Recombinant, Human Monoclonal
 Anti-GITR

Catalog #: 71295-1

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:

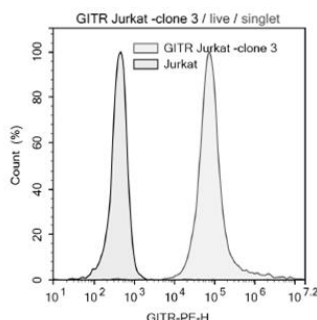
1. Cohen, A., *et al.*, *Cancer Research*.
 2006 May; **66(9)**: 4904-4912.
2. Scumpia, P.O., *et al.*, *Blood*. 2007 Oct;
110(10):3673-3681.

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
 Please visit our website at: www.bpsbioscience.com