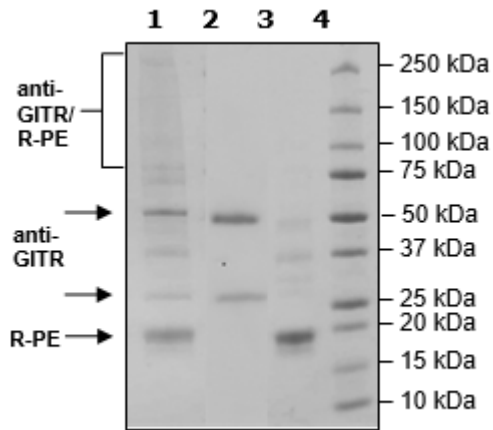


Product Information

Description:	Phycoerythrin-labeled anti-GITR recombinant human antibody recognizing the GITRL binding region of human GITR.
Label:	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, α subunit (18 kDa) and β -subunit (20 kDa), and an additional γ -subunit (34 kDa). PE is covalently attached randomly through lysines on the target protein.
Concentration:	1.68 mg/ml
Host Species:	Human
Isotype:	IgG1
Clonality:	Monoclonal
Cross Reactivity:	This antibody has not been tested for cross reactivity with any other species.
Formulated In:	8 mM Na Phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl and 20% glycerol.
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 12 months at -80°C. Avoid freeze/thaw cycles. Protect from light.
Storage:	-80°C
MW:	Heavy Chain: 51 kDa + PE, Light Chain: 26 kDa + PE
Purity:	≥90%
Purification:	Protein A affinity chromatography from HEK293 supernatants.
Assay Conditions:	For flow cytometry, we recommend incubating cells with 20 μ g/ml of anti-human GITR antibody, PE-labeled for one hour at 4°C, followed by two washes with FACS buffer.
Applications:	Useful for labeling cells expressing GITR for flow cytometry and immunofluorescence microscopy.

Quality Control Data

4-20% SDS-Page Coomassie Staining



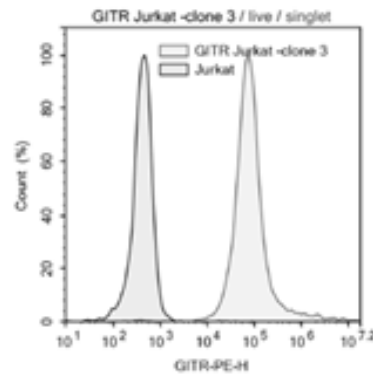
Lane 1: 8 µg anti-GITR/R-PE

Lane 2: 4 µg anti-GITR

Lane 3: 4 µg R-PE

Lane 4: Protein Marker

FACS Assay



	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.