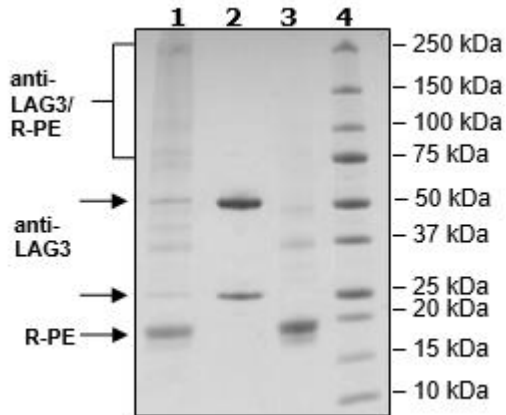


Product Information

Description:	R-Phycoerythrin-labeled recombinant human Anti-LAG3 antibody recognizing the LAG3 binding region.
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:	2.29 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 on ice in the dark for 30 minutes with 1 $\mu\text{g}/10^6$ cells of RPE anti-LAG3 antibody.
Applications:	Useful for labeling cells expressing LAG3 for flow cytometry and immunofluorescence microscopy.

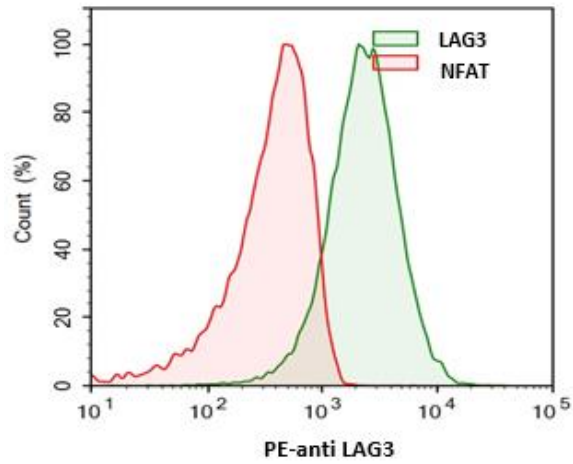
Quality Control Data

4-20% SDS-Page Coomassie Staining



Lane 1: 8 μ g anti-LAG3/R-PE
Lane 2: 4 μ g anti-LAG3
Lane 3: 4 μ g R-PE
Lane 4: Protein Marker

FACS Assay



Samples	
■	LAG3/NFAT Reporter - Jurkat
■	NFAT Reporter - Jurkat

Control Jurkat cells and LAG3-expressing stable Jurkat cells were treated with human PE anti-LAG3 antibody.