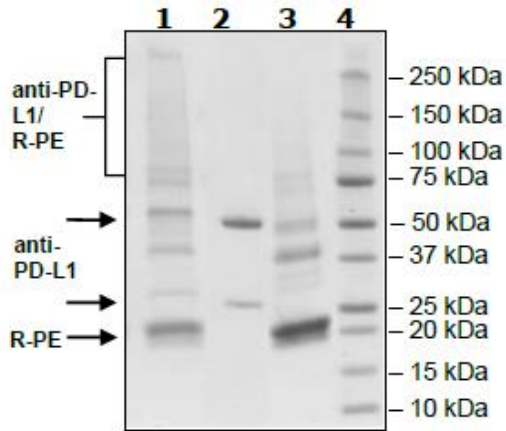


Product Information

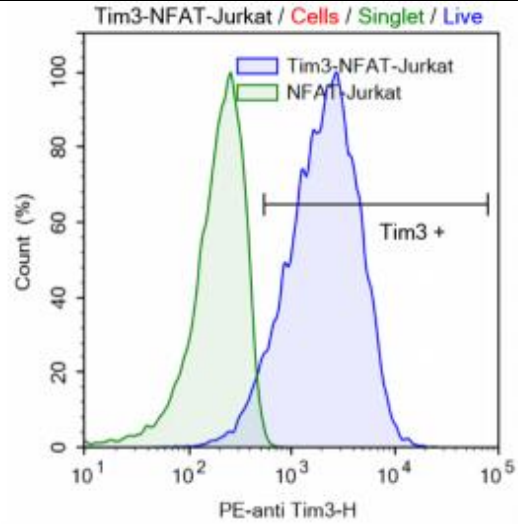
Description:	R-Phycoerythrin – labeled recombinant human Anti-TIM-3 antibody recognizing human TIM3 protein.
Label:	R-Phycoerythrin (PE) is an oligomeric protein complex (270 kDa) from red algae that exhibits intensely bright red-orange fluorescence with high quantum yields. The complex consists of six 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:	2.00 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, Light Chain: 26 kDa
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 PE-anti-TIM-3 antibody in 100 μ l of FACS buffer for 20 min at 4°C in dark.
Applications:	Useful for labeling cells expressing TIM-3 for flow cytometry and immunofluorescence microscopy discovery.

Quality Control Data

4-20% SDS-Page Coomassie Staining



FACS Assay



Samples	
■	TIM-3-NFAT-Jurkat
■	NFAT-Jurkat