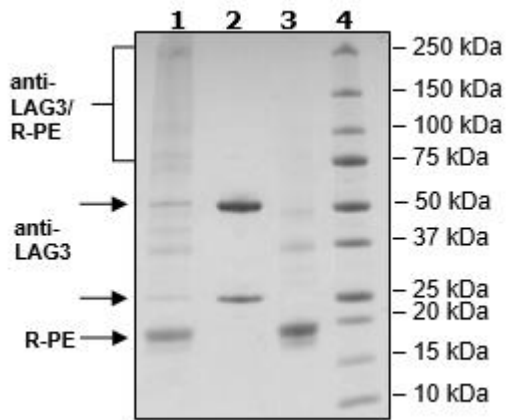


Product Information

| | |
|---------------------------|--|
| 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



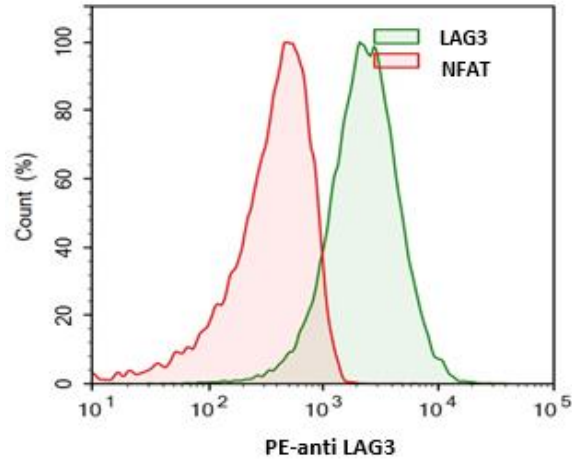
Lane 1: 8 µg anti-LAG3/R-PE

Lane 2: 4 µg anti-LAG3

Lane 3: 4 µg 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.