

Data Sheet

Anti-LAG3 Antibody, PE-Labeled

Recombinant, Human Monoclonal
Catalog #: 71226-2
Lot #: 160907 **Conc.:** 2.29 mg/ml

Host Species | Isotype:
 Human | IgG1

Formulation: 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. Avoid freeze/thaw cycles. Store
 diluted antibody with carrier protein (BSA 0.1
 – 0.5%) at 4°C, stable for 2 weeks.

References:

1. Byrnes, et al., Cytometry A. 2012
 Sep;81(2):165-8.
2. Di Carlo, et al., J. Pathol. 2005 Jan;
 205(1):82-91.

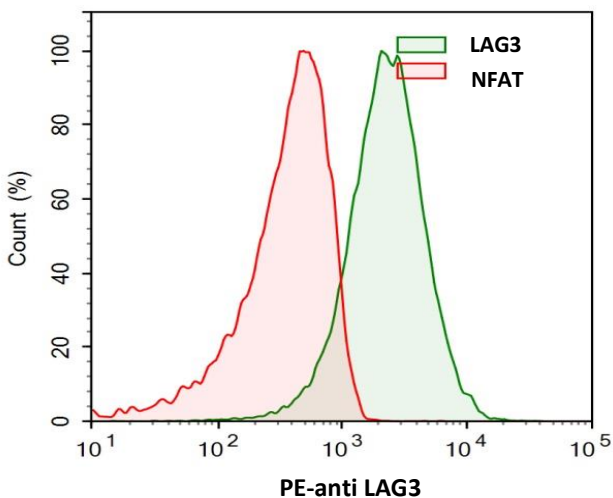
Description: R-Phycoerythrin-labeled recombinant human Anti-LAG3 antibody recognizing the LAG3 binding region.

Assay Conditions: For flow cytometry, we recommend incubating cells on ice in the dark for 30 minutes with 1 µg/10⁶ cells of RPE anti-LAG3 antibody.

Purification: Protein A affinity chromatography from HEK293 supernatants.

Application: Useful for labeling cells expressing LAG3 for flow cytometry and immunofluorescence microscopy.

Quality Assurance



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

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

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