

## Data Sheet

### Anti-LAG3 Antibody, PE-Labeled

Recombinant, Human Monoclonal  
**Catalog #:** 71226-1  
**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.

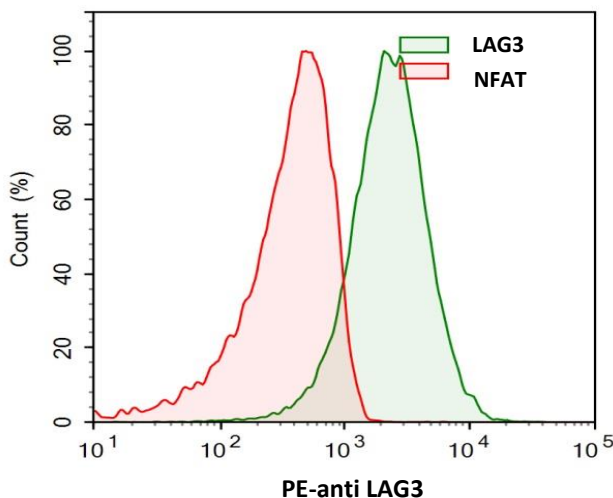
**Description:** R-Phycoerythrin-labeled  
 recombinant human Anti-LAG3 antibody  
 recognizing the LAG3 binding region.

**Assay Conditions:** For flow cytometry, we  
 recommend incubating cells on ice in the dark  
 for 30 minutes with 1 µg/10<sup>6</sup> 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](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)