

Data Sheet

Anti-LAG3, Neutralizing Antibody

Recombinant, Human Monoclonal
Catalog #: 71219
Lot #: 161012 **Conc.:** 8.79 mg/ml

Host Species | Isotype:
 Human | IgG1

Formulation: 8 mM 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: Neutralizing recombinant human antibody recognizing the Lag-3 binding region.

EC50: 0.58 µg/ml BPS LAG3 Reporter Cell Line (BPS Cat. #71278).

Assay Conditions: Anti-LAG3 Neutralizing Antibody was added at various concentrations to the LAG-3/NFAT Reporter Jurkat cells (BPS Cat. #71278) co-cultured with superantigen-loaded Raji cells and luciferase activity was measured using ONE-Step Luciferase Assay System (BPS Cat. #60690)

Purification: Protein A affinity chromatography from HEK293 supernatants.

Application: Useful as a reference antibody for screening LAG-3 neutralizing antibodies.

Quality Assurance

Neutralizing Anti-LAG3 Activity

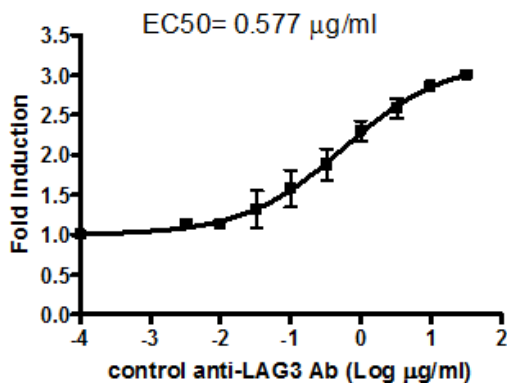


Figure 1. Dose response of anti-LAG3 neutralizing antibody in LAG-3/NFAT Reporter-Jurkat cells (BPS Cat #71278).

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