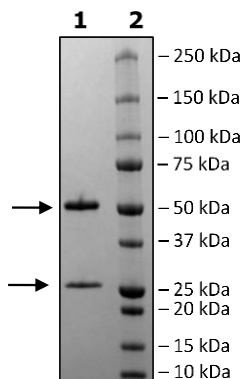


Product Information

Description:	Neutralizing recombinant human antibody recognizing the Lag-3 binding region.
Concentration:	2.09 mg/ml
Species:	Human
Formulated In:	8 mM phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
Expression System:	HEK293
Isotype:	IgG1
Clonality:	Monoclonal
Format:	Aqueous buffer solution
Stability:	At least 12 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
MW:	Heavy Chain: 50 kDa + glycans; Light Chain: 23 kDa
Purity:	≥90%
Purification:	Protein A affinity chromatography from HEK293 supernatants.
Assay Conditions:	Assay was performed according to the LAG3:FGL1[Biotinylated] Inhibitor Screening Assay Kit (BPS Bioscience #79789) with Anti-LAG3 titrated from 500 nM to 0 nM
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



LAG3:FGL1[Biotinylated] Activity

