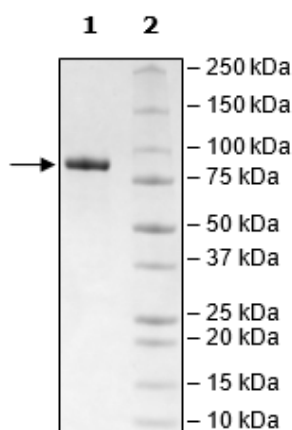


Product Information

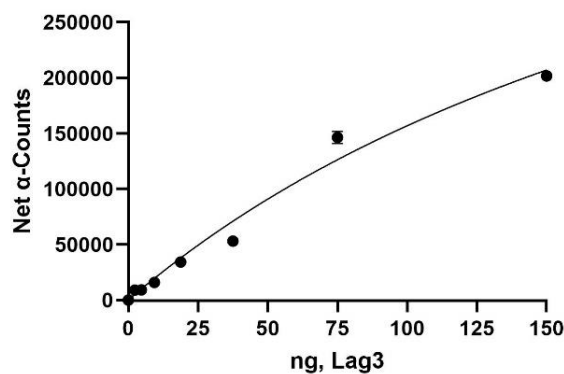
Construct:	LAG3 (23-450-Fc(IgG1)-Avi)-(Biotin)
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.
Concentration:	1.36 mg/ml
Species:	Human
Formulated In:	20 mM NaOAc, pH 5.5, 500 mM NaCl, 0.02% Tween 20 and 10% glycerol
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_002286
MW:	75 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	≥90%
Aggregation:	<10%
Assay Conditions:	Assay was conducted in 1x Immuno Buffer 1 (BPS Bioscience #79311). 5 µl FGL1 His from Cloud Clone Corp. #RPD022Hu01 P20211110840 was used with a final concentration of 100 nM along with 5 µl of titrated LAG3 (CD223), Biotin-labeled (Human) HiP™ (BPS Bioscience #71147) titrated from 150-0 ng/rxn and incubated at room temperature for 1 hour. 10 µl 1:250-fold diluted Nickel Chelate acceptor beads Perkin Elmer #AL108C were added to the plate and incubated at room temperature for 30 min. 10 µl 1:125 fold diluted Strep Donor Beads Perkin Elmer #6760002S were added and incubated at room temperature for 1 hour. A-counts were read using a Tecan infinite M1000 Pro plate reader.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

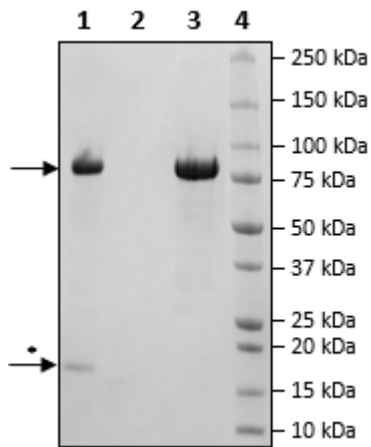
4-20% SDS-Page Coomassie Staining



LAG3-FGL1 Activity



Biotin-Avidin Pulldown



1. Beads
2. Flow thru
3. Control
4. Standards

* Avidin from beads.

Gel Filtration Trace

