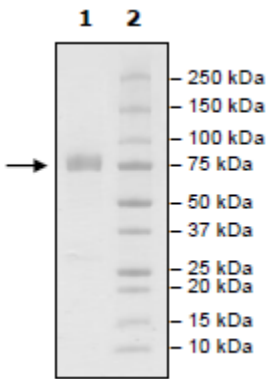


Product Information

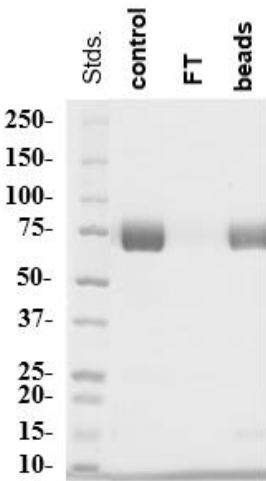
Construct:	PD-L1(Mouse)(19-238-mIgG2aFc-Avi-Biotin)
Label:	This protein was biotinylated using Avi-Tag™ technology. Biotinylation confirmed to be >90%.
Concentration:	1.32 mg/ml
Species:	Mouse
Formulated In:	8 mM Phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol.
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_021893
MW:	54 kDa
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Aggregation:	<10% as confirmed by gel filtration.
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining



Biotin-Avidin Pulldown



Gel Filtration Trace

