

PD-L1, Fc Fusion, PE-Labeled (Mouse) Recombinant

Catalog: 71268
Lot: 230420

Product Information

Description:	Mouse PD-L1 (Programmed Death-Ligand 1), also known as CD274 or B7-H1, encompassing amino acids 18 to 238. This construct has a mouse IgG2a Fc domain on the C-terminus and has been labeled with PE. The recombinant protein was affinity purified.
Species:	Mouse
Construct:	mPD-L1 (18-238-Fc(mIgG2a))-(PE)
Concentration:	1.55 mg/ml
Expression System:	HEK293
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	51 kDa + glycans + PE
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	NM_021893
Label:	R-Phycoerythrin (PE) is an oligomeric protein complex (270 kDa) from red algae that exhibits intensely bright red-orange fluorescence with high quantum yields. The complex consists of six heterodimers, α subunit (18 kDa) and β -subunit (20 kDa), and an additional γ -subunit (34 kDa). PE is covalently attached randomly through lysines on the target protein.
Stability:	At least 6 months at -80°C. Protect from light.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

