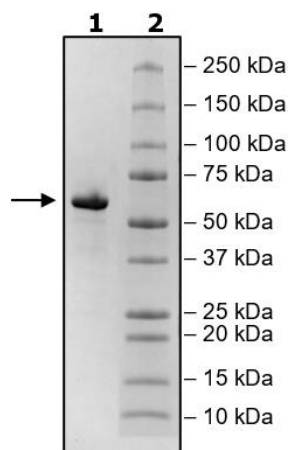


Product Information

Description:	Recombinant mouse CD73 (5'-nucleotidase), encompassing amino acids 29-549. This construct contains a C-terminal Avi-His-tag (6xHis). This recombinant protein was affinity purified.
Species	Mouse
Construct:	CD73 (Mouse) (29-549-Avi-His)
Concentration:	1.12 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:	61 kDa + glycans
Glycosylation	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	NM_011851
Stability:	At least 6 months at -80°C.
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.
Specific Activity:	≥20 nmol/min/μg
Assay Conditions:	Assay was performed according to CD73 Inhibitor Screening Assay Kit (BPS Bioscience #72055) with CD73 enzyme titrated at various concentrations.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



CD73 (Mouse) Activity

