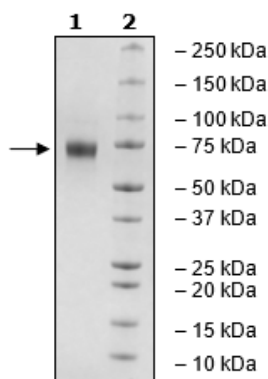


## Product Information

<b>Construct:</b>	CD39 (38-478-His)
<b>Concentration:</b>	0.50 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
<b>Expression System:</b>	HEK293
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_001776
<b>MW:</b>	52 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Purity:</b>	≥90%
<b>Specific Activity:</b>	≥3000 pmol/min/μg
<b>Assay Conditions:</b>	Assay was done according to CD39 Inhibitor Screening Assay Kit (BPS Bioscience #79278) with various amounts of CD39.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-PAGE Coomassie Staining



CD39 Activity

