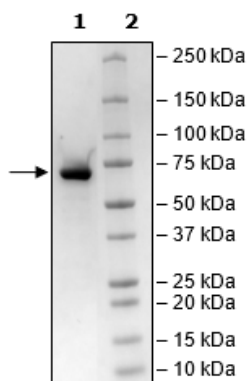


Product Information

Construct:	SARM1 (FLAG-28-724(end))
Concentration:	0.78 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 3 mM DTT, and 100 µg/ml FLAG peptide
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_015077
MW:	77 kDa
Purity:	≥90%
Assay Conditions:	Assay was done according to the SARM1 Fluorogenic Assay Kit (Hydrolase Activity) (BPS Bioscience #78217).
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



SARM1 Activity

