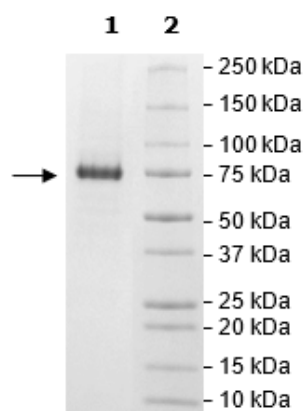


Product Information

Construct:	Sphingosine Kinase 2 (Transcript Variant 1) (His-1-654(end))
Concentration:	0.91 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% Glycerol and 250 mM imidazole
Expression System:	Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_020126
MW:	73 kDa
Purity:	≥90%
Specific Activity:	15 pmol/min/μg
Assay Conditions:	Assay was done according to the SPHK1 Assay Kit (BPS Bioscience #78026) with various amounts of SPHK2.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

