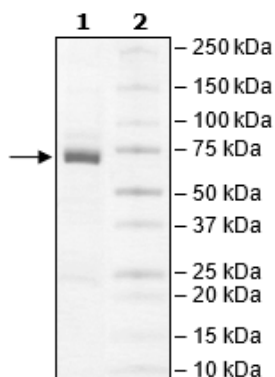


Product Information

Construct:	Sphingosine Kinase 2 (His-1-618(end))
Concentration:	0.29 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT, variable amounts of Imidazole, and 20% Glycerol.
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_001204160
MW:	70 kDa
Purity:	≥90%
Specific Activity:	20 pmol/min/μg
Assay Conditions:	The activity of SPHK2 was measured with increasing amounts of protein according to the protocol described in SPHK1 Assay Kit (BPS Bioscience #78026).
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



SPHK2 Activity

