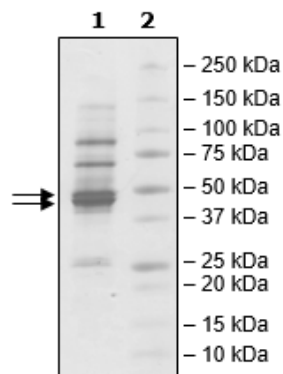


## Product Information

<b>Construct:</b>	Sphingosine-1-phosphate lyase1 (59-568(end)-His)
<b>Concentration:</b>	0.37 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	50 mM Tris, pH 8.0, 500 mM NaCl, 100 mM imidazole, 3 mM methionine, 1 mM pyridoxal 5'-phosphate monohydrate, 1 mM TCEP, and 10% glycerol
<b>Expression System:</b>	Sf9
<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_003901
<b>MW:</b>	58 kDa
<b>Purity:</b>	54%
<b>Assay Conditions:</b>	S1PL-1 reaction was done in 100 mM Phosphate buffer (pH=7.2) containing 200 μM Fluorogenic substrate, 250 μM PLP and S1PL-1 for 2 hours at 37°C. Fluorescence signal was read at Ex360/Em450.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-Page Coomassie Staining



S1PL-1 Activity

