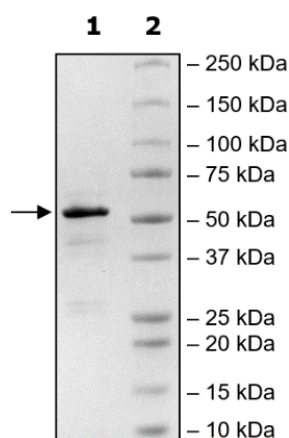


Product Information

Construct:	Sirtuin 3 (GST-102-399(end))
Concentration:	0.33 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution.
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_012239
MW:	59 kDa
Purity:	89%
Specific Activity:	3.7 pmol/min/μg
Assay Conditions:	25 mM Tris/HCl, pH 8.0, 137 mM NaCl, 2.7 mM KCl, 1 mM MgCl ₂ , and 0.1 mg/ml BSA, 20 μM Fluorogenic HDAC/SIRT substrate 1 (BPS Bioscience #50032), 500 μM NAD ⁺ and Sirtuin 3 (31.25 ng - 2000 ng). Incubation condition: 1 hour at 37°C followed by treatment with 50 μl of SIRT developer (BPS Bioscience #50089) for 15 min at room temperature. Fluorescence intensity is measured at ex360/em460.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



SIRT3 Activity

