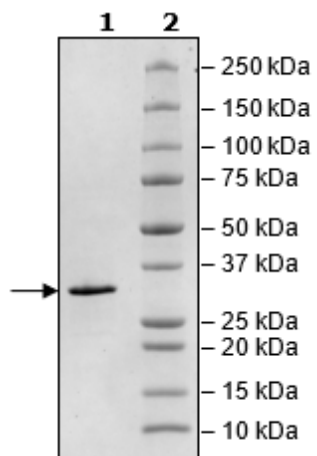


Product Information

Description:	Recombinant human Sirtuin 2 (SIRT2), encompassing amino acids 50-356. This construct contains a C-terminal His tag (6xHis). This recombinant protein was affinity purified.
Species	Human
Construct:	Sirtuin 2 (50-356-His)
Concentration:	2.29 mg/ml
Expression System:	<i>E. coli</i>
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	36 kDa
Genbank Accession:	NM_012237
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Specific Activity:	1 pmol/min/μg
Assay Conditions:	Sirtuin 2 activity was measured using the SIRT2 (Sirtuin2) Fluorogenic Assay Kit (BPS Bioscience #50087) with SIRT2 titrated at various concentrations.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



SIRT2 Activity

