



11526 Sorrento Valley Rd SteA2
San Diego Ca 92121
Tel: 1.858.202.1401
1.858.829.3082
Fax: 1.858.259.5734
Email: info@bpsbioscience.com

Data Sheet

Sirtuin1

human, recombinant, N-terminal GST tag
Catalog #: 50012

Formulated in:

25 mM Tris-HCl, pH 8.0, 100 mM NaCl, 0.05% Tween-20 and 20% glycerol.

Stability:

>6 months at -80°C

References:

1. Kolthur-Seetharam, U. et al (2006) Cell Cycle 5, 873-877.
2. Solomon, J.M. et al (2006) Mol. Cell. Biol. 26, 28-38.

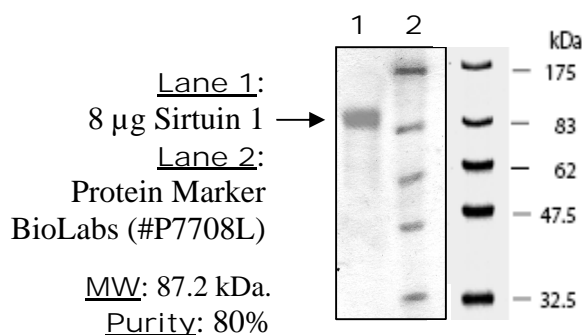
Description: Human Sirtuin 1 (GenBank Accession No. NM_012238), amino acids 193- 741 with N-terminal GST tag, MW= 87.2 kDa, expressed in E.coli expression system.

Specific Activity: 8 pmol/min / μ g. Assay condition: 25 mM Tris/Cl, pH8.0, 137 mM NaCl, 2.7 mM KCl, 1 mM MgCl₂, and 0.1 mg/ml BSA, 30 μ M Biomol substrate (Catalog number KI-177), 200 μ M NAD⁺ and 20 ng/ μ l Sirtuin 1. Incubation condition: 30 min at 30°C.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

10% SDS-PAGE Coomassie staining



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** or **1.858.829.3082**, Fax **1.858.259.5734**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com