

Data Sheet

Sirtuin3, GST-Tag

Human, Recombinant, N-terminal GST tag

Catalog #: 50014

Lot #: 181005

Conc.: 0.23 mg/ml

Formulated in: 40mM Tris-HCl pH8.0, 110mM NaCl, 2.2mM KCl, and 20% glycerol

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

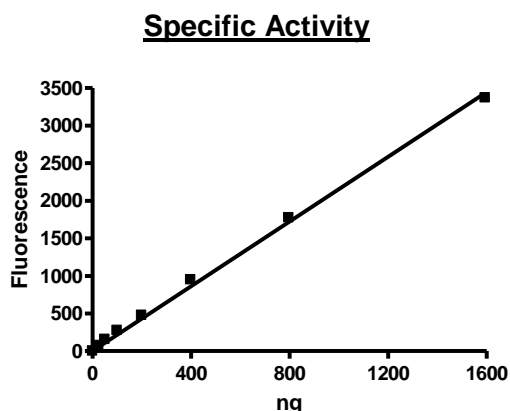
1. Scher, M.B. *et al.*, *Genes Dev.* **21 (8)**, 920-928 (2007)
2. Bellizzi, D. *et al.*, *Genomics* **89 (1)**, 143-150 (2007)

Description: Human Sirtuin 3 (GenBank Accession No. NM_012239), amino acids 102-399 with N-terminal GST tag, MW= 59 kDa, expressed in *E. coli* expression system.

Specific Activity: 3.7 pmole/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 BPS HDAC substrate 1 (Catalog #50032), 500 μM NAD^+ and Sirtuin 3 (16 – 2000 ng). Incubation condition: 30 min at 37°C followed by developing for 15 min at room temperature.

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

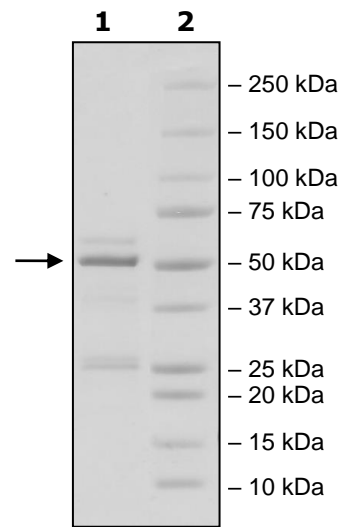
Quality Assurance



**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
3 μg SIRT3
Lane 2:
Protein Marker

MW: 59kDa
Purity: 63%



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

To place your order, please contact us by Phone **1.858.829.3082**, Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com