

Data Sheet

Sirtuin 5

Human, recombinant, N-terminal GST tag

Catalog #: 50016

Lot#: 130110 **Conc.:** 0.94 mg/ml

Formulated in: 45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 18 mM glutathione and 10% glycerol.

Stability: >6 months at -80°C

References:

1. Mahlknecht, U. *et al.*, (2006) *Cytogenet. Genome Res.* **112**, 208-212.
2. Frye, R.A. (2000) *Biochem. Biophys. Res. Commun.* **273**, 793-798.

Description:

Human Sirtuin 5 (GenBank Accession No. NM_012241), full length with N-terminal GST tag, MW= 59.8 kDa, expressed in *E. coli* expression system.

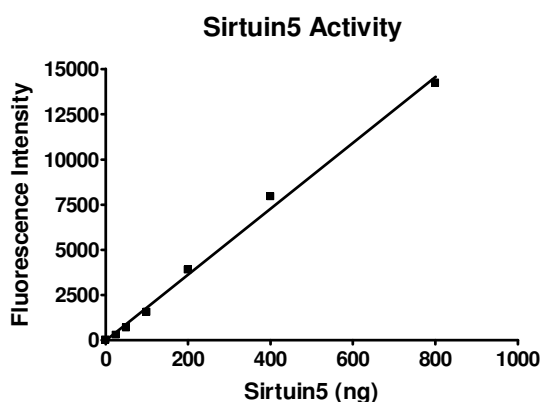
Specific Activity: 3.6 pmole/min/μg

Assay Conditions: 50 mM Tris-HCl, pH 8.0, 137 mM NaCl, 2.7 mM KCl, 1 mM MgCl₂, 0.05% Tween 20, 500 μM NAD⁺, 200 μM Sirtuin 5 substrate, and Sirtuin 5. Incubate for 30 minutes at 37°C. Fluorescence intensity is measured at exc360/em460.

Application:

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

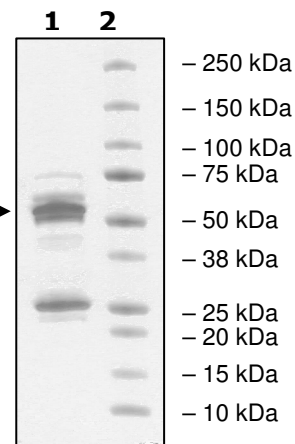
Quality Assurance



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

Lane 1:
 6 μg Sirtuin 5
Lane 2:
 Protein Marker

MW: 59.8 kDa
Purity: ≥50%



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