



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](mailto:info@bpsbioscience.com)

## Data Sheet

### Sirtuin4

human, recombinant, N-terminal GST tag  
Catalog #: 50015

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.Frye, R.A. (2000) Biochem. Biophys. Res. Commun. 273, 793-798.

Description:

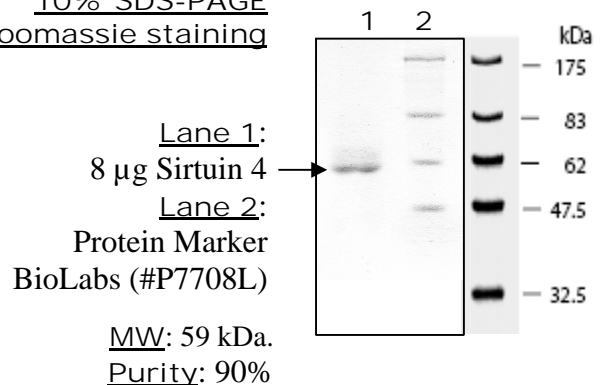
Human Sirtuin 4 (GenBank Accession No. NM\_012240), amino acids 25-314 with N-terminal GST tag, MW= 59 kDa, expressed in E.coli expression system.

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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)