



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

### **Sirtuin 5**

Human, recombinant, N-terminal GST tag  
Catalog #: 50016

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

**Description:**

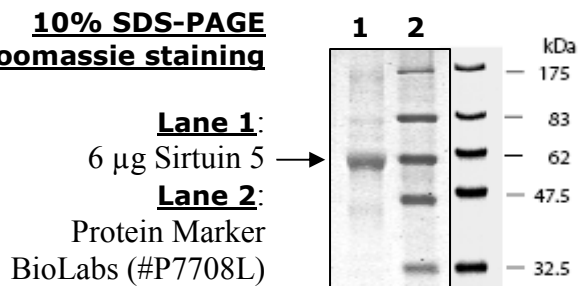
Human Sirtuin 5 (GenBank Accession No. NM\_012241), full length with N-terminal GST tag, MW= 59.8 kDa, expressed in E.coli expression system.

**Application:**

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

### **Quality Assurance**

**10% SDS-PAGE  
Coomassie staining**



**Lane 1:**

6 µg Sirtuin 5

**Lane 2:**

Protein Marker

BioLabs (#P7708L)

**MW:** 59.8 kDa.

**Purity:** >75%

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)