



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

Sirtuin 7

Human, recombinant, C-terminal Flag tag

Catalog #: 50018

Lot# 130111-GC1 **Conc.:** 0.13 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 0.04% Tween-
20, and 20% glycerol.

Stability: >6 months at -80°C

References:

1. Grob A. *et al.*, *J. Cell. Sci.* **122** (PT 4), 489-498 (2009).
2. Ford E. *et al.*, *Genes Dev.* **20** (9), 1075-1080 (2006).

Description:

Human Sirtuin 7 (GenBank Accession No. NM_016538), full length, MW= 46 kDa, with C-terminal Flag tag expressed in an *E. coli* cell expression system.

Application:

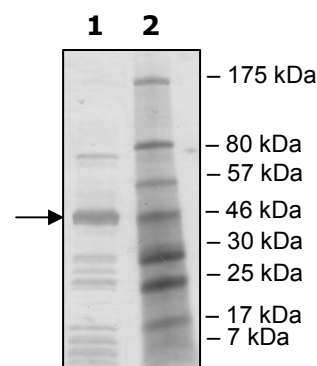
Useful for screening inhibitors, and selectivity profiling.

Quality Assurance

4-20% SDS-PAGE
Coomassie staining

Lane 1:
3 µg Sirtuin 7
Lane 2:
Protein Marker

MW: 46 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