

Data Sheet

SMYD3

Human, recombinant, N-terminal GST tag

Catalog #: 51015

Lot#: 130213 **Conc.:** 0.13 mg/ml

Formulated in: 40 mM Tris, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 16 mM glutathione, and 20% glycerol.

Stability: >6 months at -80°C

References:

1. Wang S.Z. *et al.*, *BMB Rep* **41 (4)**, 294-299 (2008).
2. Silva F.P. *et al.*, *Oncogene* **27 (19)**, 2686-2692 (2008).

Description:

Human SMYD3 (GenBank Accession No. NM_022743), amino-acids 35-end, with N-terminal GST tag, MW=64 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

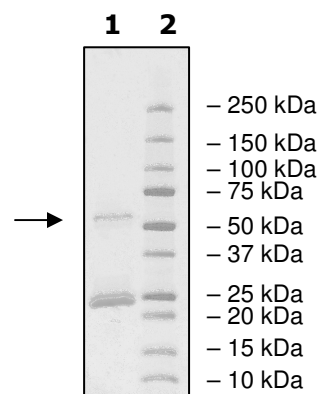
Application:

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

Quality Assurance

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

Lane 1:
2 µg SMYD3
Lane 2:
Protein Marker
MW: 64 kDa
Purity: ≥26%



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