

Data Sheet

SMYD3

Human, recombinant, N-terminal GST tag

Catalog #: 51015

Lot#: 110504-20 Conc.: 0.06 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 16 mM
Glutathione, 20% Glycerol and 3mM DTT.

Stability: >6 months at -80°C

References:

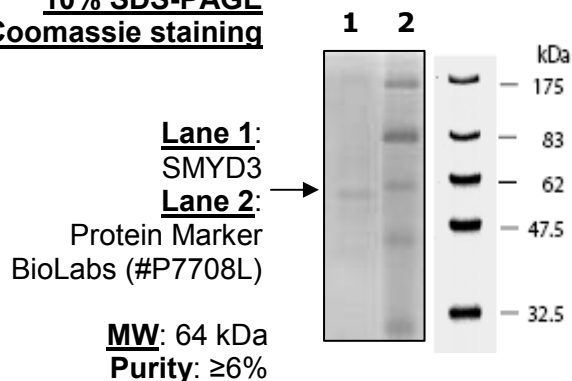
1. Wang S.Z. et al., BMB Rep **41 (4)**, 294-299 (2008).
2. Silva F.P. et al., Oncogene **27 (19)**, 2686-2692 (2008).

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

**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 **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com