

## Data Sheet

### SMYD3

Human, recombinant, N-terminal GST tag

**Catalog #:** 51015

Lot#: 110504-5      Conc.: 0.06 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0,  
110 mM NaCl, 2.2 mM KCl, 4 mM  
Glutathione, 20% Glycerol and 3mM DTT.

**Stability:** >6 months at -80°C

#### References:

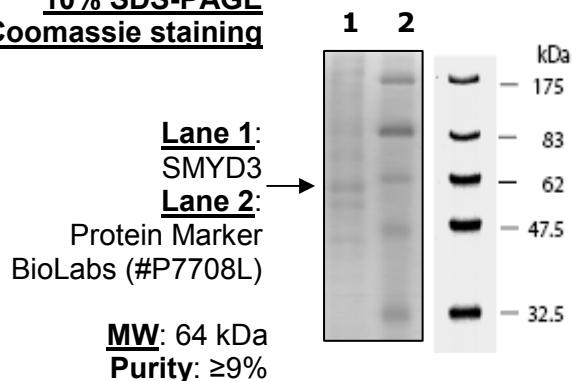
1. Wang S.Z. et al., BMB Rep **41 (4)**, 294-299 (2008).
2. Silva F.P. et al., Oncogene **27 (19)**, 2686-2692 (2008).

**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](mailto:info@bpsbioscience.com)

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