

Data Sheet

SMYD2, FLAG-Tag

Human, Recombinant, N-terminal FLAG-tag

Catalog #: 51014

Lot#: 180622 **Conc.:** 0.60 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween20, 20% glycerol, 0.2mM TCEP, 80 µg/ml FLAG peptide.

Stability: At least 6 months at -70°C. Avoid freeze/thaw cycles.

References:

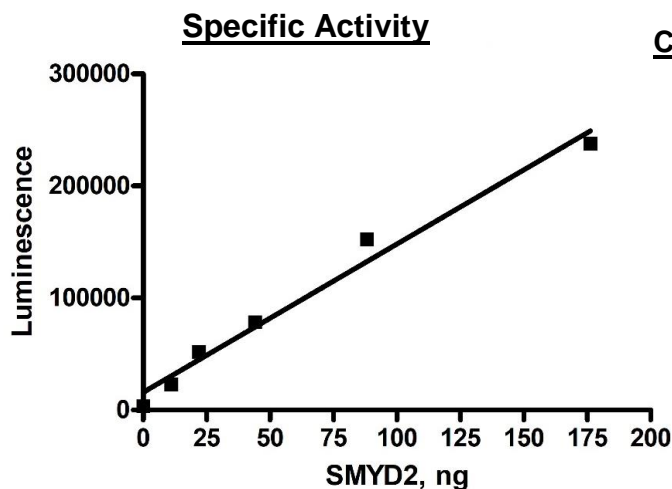
1. Abu-Farha M. et al., Mol. Cell Proteomics **7 (3)**, 560-572 (2008).
2. Brown M.A. et al., Mol. Cancer **5 (26)**, (2006).

Description: Human SMYD2, also known as KMT3C, SET And MYND Domain Containing 2 and Lysine N-Methyltransferase 3C, GenBank Accession No. NM_020197, a.a. 2-433(end) with N-terminal FLAG-tag, expressed in a Baculovirus infected Sf9 cell expression system, MW= 51 kDa.

Specific Activity: .004 pmole/min/µg
 Assay Conditions: For standard, H3, 21-44 K36 (1Me)-Biotin was coated onto a neutravidin plate at varying concentrations. For test lots, H3,21-44-Biotin was coated onto a neutravidin plate at 0.4 µg/ml. Reaction mixture contained 50 mM Tris buffer pH 8.8, 5 mM MgCl₂, 1 mM TCEP, 0.05% Tween-20, 10 µM S-adenosylmethionine, and various amounts of SMYD2—incubated for overnight at room temperature. Followed by 1 hour incubation with primary antibody against H3 K36 (1Me), and 30 minute incubation with secondary antibody. Luminescence was developed with HRP substrate A and B (Pierce).

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

Quality Assurance



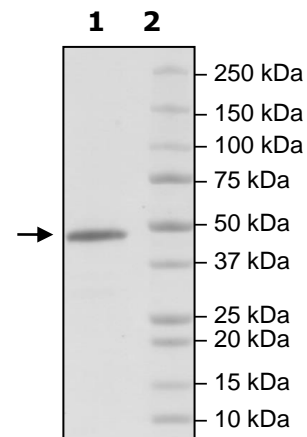
4-20% SDS-PAGE
Coomassie staining

Lane 1:
SMYD2

Lane 2:
Protein Marker

Purity: ≥90%

MW: 51 kDa



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** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com