

## Data Sheet

### PRMT2, FLAG-Tag

Human, Recombinant, N-terminal FLAG-Tag,  
**Catalog #:** 100471  
**Lot #:** 190827-2      **Conc.:** 0.36 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween20, 100 µg/ml FLAG peptide, 3mM DTT, 20% glycerol.

**Stability:** At least 6 months at -80°C. Avoid freeze/thaw cycles.

#### References:

1. Blythe, S.A., *et al.*, *Developmental Cell*. 2010 Aug; **19(2)**:220-231.
2. Yoshimoto, T., *et al.*, *Experimental Cell Research*. 2006 Jul; **312(11)**: 2040-2053.

**Description:** Human Protein Arginine N-Methyltransferase 2 (PRMT2), also known as Histone-arginine N-methyltransferase PRMT2; HMT1, and HRMT1L1, GenBank Accession No. NM\_206962, a.a. 2-433(end), with an N-terminal FLAG-tag, expressed in an SF9 insect cell expression system. MW = 50 kDa.

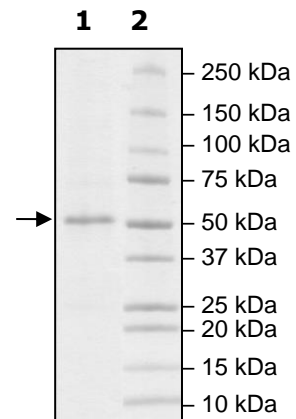
## Quality Assurance

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

**Lane 1:**  
PRMT2  
**Lane 2:**  
Protein Marker

**Purity:** ≥90%

**MW:** 50 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](mailto:info@bpsbioscience.com)

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