

Data Sheet

JMJD6

Human, recombinant, N-terminal GST-tag

Catalog #: 50140

Lot#: 121218-GC **Conc.:** 1.87 mg/ml

Formulated in: 40 mM Tris-HCl, pH 7.5, 110 mM NaCl, 2.2 mM KCl, 0.05% Tween-20, and 20% Glycerol.

Stability: >6 months at -80 °C

References:

1. Hahn, P., *et al.* (2010). *PLoS One* **5(10):**e13769.
2. Webby, C.J., *et al.* (2009). *Science* **325(5936):**90-3.
3. Chang, B., *et al.* (2007). *Science* **318(5849):**444-7.

Description:

Human JMJD6 (Jumonji domain containing 6), also known as PTDSR (GenBank Accession No. NM_001081461), amino-acids 2-414 (end) with an N-terminal GST-tag, MW = 74 kDa, expressed in a Baculovirus-infected Sf9 cell expression system.

Application:

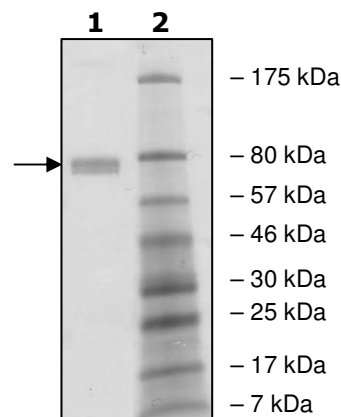
Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
2 µg JMJD6
Lane 2:
Protein Marker

MW: 74 kDa
Purity: ≥85%



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