

## Data Sheet

### JMJD2B (KDM4B), FLAG-tag

Human, Recombinant, C-terminal FLAG-tag

**Catalog #:** 50104

**Lot #:** 170309-1      **Conc.:** 0.16 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 80 µg/mL FLAG peptide and 20% glycerol

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

**References:**

1. Fodor, B.D., *et al.* (2006). *Genes Dev.* **20(12)**: 1557–1562.
2. Beyer S., *et al.* (2008). *J. Biol. Chem.* **283(52)**:36542-52.

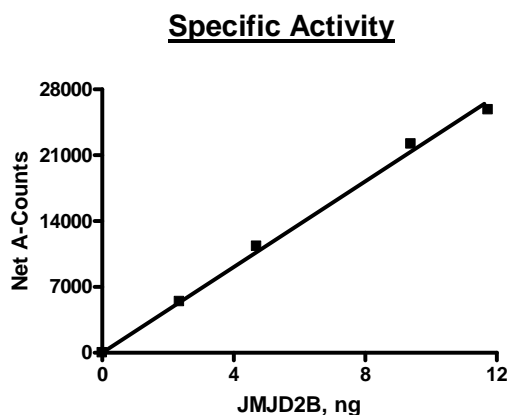
**Description:** Human JMJD2B, also known as JHDM3B and KDM4B (GenBank Accession No. NM\_015015) (amino-acids 1-500) with C-terminal FLAG tag, MW=57 kDa, expressed in Sf9 cells using a Baculovirus expression system.

**Specific Activity:**  $\geq 0.23$  pmol/min/µg.

Assay conditions: 10µl reaction mix containing assay buffer with 20 mM HEPES (pH 7.4), 50 mM NaCl, 1mM TCEP, 500 µM  $\alpha$ -ketoglutarate, 25 µM iron, 2mM ascorbic acid, 0.01% Tween-20, 0.5 µM biotinylated peptide substrate, and JMJD2B (2 – 10ng) added to the wells. Add antibody against dimethylated K<sub>9</sub> peptide. Incubate for 30 min, then Streptavidin-conjugated secondary antibody followed by Alpha Screening detection.

**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 JMJD2B  
**Lane 2:**  
Protein Marker

**MW:** 57 kDa  
**Purity:**  $\geq 61\%$

