

---

## JMJD2B Protein

(Catalog No. E24023)

---

### Description

Human JMJD2B, also known as JHDM3B and KDM4B (GenBank Accession No. NM\_015015), a.a. 2-500 with N-terminal GST-tag.

### Molecular Weight (kDa)

83 kDa

### Formulation

40 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 20% glycerol, 16 mM glutathione, and 3mM DTT.

### Source

Expressed in Sf9 cells using a Baculovirus expression system.

### Size

20 µg

### Storage Temp & Conditions

>6 months at -80°C

### Purification Notes

Purity >40%

### Application Notes

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

### Ordering Information:

Products	Size	Cat. No.
JMJD2B Protein	20 µg	E24023 -1

*This product is for research purposes only. Not intended for use in diagnostic procedures.*