
JMJD3 Protein

(Catalog No. E24026)

Description

Human JMJD3/KDM6B histone demethylase (a.a. 1043-end), Genbank Accession No. NM_001080424, with C-terminal DDDK FLAG-tag.

Molecular Weight (kDa)

71 kDa

Formulation

45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 90 µg/ml Flag peptide, 10% glycerol, and 3 mM DTT.

Source

Expressed in Sf9 cells via a baculovirus expression system

Size

20 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Purity>50%

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
JMJD3 Protein	20 µg	E24026 -1

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