
JMJD2A Protein

(Catalog No. E24022)

Description

Human JMJD2A, also known as JHDM3A and KDM4A (GenBank Accession No. NM_014663) (amino-acids 1-350) with N-terminal HIS-tag.

Molecular Weight (kDa)

42 kDa

Formulation

40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 250 mM imidazole, and 20% Glycerol.

Source

Expressed in an E.coli expression system

Size

20 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Purity >55%

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
JMJD2A Protein	20 µg	E24022 -1

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