
JMJD2D Protein

(Catalog No. E24025)

Description

Human JMJD2D (GenBank Accession No. NM_018039) (amino-acids 2- 354) with N-terminal GST tag.

Molecular Weight (kDa)

67 kDa

Formulation

40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 4 mM Glutathione, 20% Glycerol and 3 mM DTT.

Source

Expressed in a Baculovirus infected Sf9 cell expression system.

Size

20 µg

Storage Temp & Conditions

>6 months at -80°C

Purification Notes

Purity >30%

Application Notes

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

Ordering Information:

Products	Size	Cat. No.
JMJD2D Protein	20 µg	E24025 -1

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