

---

## JMJD2C Protein

(Catalog No. E24024)

---

### Description

Human JMJD2C (GenBank Accession No. BC143571) (amino-acids 2-372) with N-terminal GST tag, MW=69 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

### Molecular Weight (kDa)

69 kDa

### Formulation

45 mM Tris-HCl, pH8.0, 124 mM NaCl, 2.4 mM KCl, 18 mM glutathione, 10% 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 >80%

### Application Notes

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

### Ordering Information:

| Products       | Size  | Cat. No.  |
|----------------|-------|-----------|
| JMJD2C Protein | 20 µg | E24024 -1 |

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