

---

## JARID1C Protein

(Catalog No. E24021)

---

### Description

Human JARID1C, also known as SMCX (Smcy homolog, X-linked), KDM5C, or XE169, GenBank Accession No. NM\_004187), full length [a.a. 2-1560(end)] with N-terminal FLAG- and His-tags.

### Molecular Weight (kDa)

181 kDa

### Formulation

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

### Source

Expressed in Sf9 cells via 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.
JARID1C Protein	20 µg	E24021 -1

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