

---

## SMYD3 Protein

(Catalog No. E24036)

---

### Description

Human SMYD3 (GenBank Accession No. BC031010), full length, with N- terminal FLAG-tag.

### Molecular Weight (kDa)

50 kDa

### Formulation

40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 80 µg/ml FLAG peptide, 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>90%

### Application Notes

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

### Ordering Information:

Products	Size	Cat. No.
SMYD3 Protein	20 µg	E24036 -1

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