

---

## EZH2 Protein

(Catalog No. E24031)

---

### Description

Complex of human EZH2 (GenBank Accession No NM\_004456), (amino acids 2-end) with N-terminal His-GST tag.

### Molecular Weight (kDa)

114 kDa

### Formulation

40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT, 16 mM glutathione and 20% Glycerol.

### Source

Expressed in a Sf9 baculovirus expression system

### Size

20 µg

### Storage Temp & Conditions

>6 months at -80°C

### Purification Notes

Purity>65%

### Application Notes

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

### Ordering Information:

Products	Size	Cat. No.
EZH2 Protein	20 µg	E24031 -1

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