

EZH1 Protein

Recombinant full-length human protein expressed in Sf9 cells

Catalog # E395-30G

Lot # Q2465-7

Product Description

Recombinant full-length human EZH1 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The EZH1 gene accession number is [BC015882](#).

Gene Aliases

KIAA0388

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, and 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

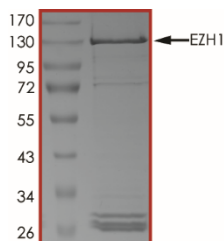
Scientific Background

EZH1 (enhancer of zeste homolog 1) belongs to the polycomb group (PcG) protein and acts as the catalytic subunit of the PRC2/EED-EZH1 complex, which methylates 'Lys-27' of histone H3, leading to transcriptional repression of target genes. EZH1 is ubiquitously expressed. Compared to PRC2-EZH2, PRC2-EZH1 exhibits low levels of histone methyltransferase (HKMT) activity. Required for embryonic stem cell derivation and self-renewal, suggesting that it is involved in safeguarding embryonic stem cell identity.

References

1. Margueron R, et al: Ezh1 and Ezh2 maintain repressive chromatin through different mechanisms. *Mol Cell*. 32(4):503-18, 2008.
2. Shen X, et al: EZH1 mediates methylation on histone H3 lysine 27 and complements EZH2 in maintaining stem cell identity and executing pluripotency. *Mol Cell*. 32(4):491-502, 2008.

Purity



The purity of EZH1 was determined to be **>70%** by densitometry, approx. MW **130 kDa**.

EZH1 Protein

Recombinant full-length human protein expressed in Sf9 cells

Catalog #	E395-30G
Lot #	Q2465-7
Purity	>70%
Concentration	0.1 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.