

EZH2 (Y641N) Protein

Recombinant full-length human protein expressed in Sf9 cells

Catalog # E396-32EG

Lot # B2085-2

Product Description

Recombinant full-length human EZH2 (Y641N) protein was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The EZH2 gene accession number is [BC010858](#).

Gene Aliases

ENX-1, EZH1, MGC9169

Formulation

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

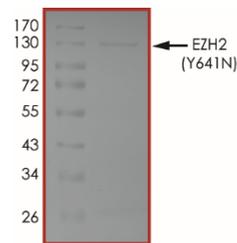
Scientific Background

EZH2 (enhancer of zeste homolog 2) belongs to the polycomb group (PcG) protein and acts as the catalytic subunit of the PRC2/EED-EZH2 complex, which methylates 'Lys-27' of histone H3, leading to transcriptional repression of target genes. EZH2 expression is associated with proliferating tissues and shares an overlapping set of target genes with EZH1. Compared to PRC2-EZH1, PRC2-EZH2 exhibits higher levels of histone methyltransferase (HKMT) activity.

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 EZH2 (Y641N) was determined to be **>80%** by densitometry, approx. MW **130 kDa**.

EZH2 (Y641N) Protein

Recombinant full-length human protein expressed in Sf9 cells

Catalog #	E396-32EG
Lot #	B2085-2
Purity	>80%
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.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.