

Recombinant SUV420H2 protein

Catalog No: 31361

Expressed In: *E. coli*

Quantity: 50 µg

Concentration: 0.07 µg/µl

Source: Human

Buffer Contents: 50 µg recombinant SUV420H2 supplied at a concentration of 0.07 µg/µl in a buffer of 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 16 mM Glutathione, 20% glycerol, and 3 mM DTT.

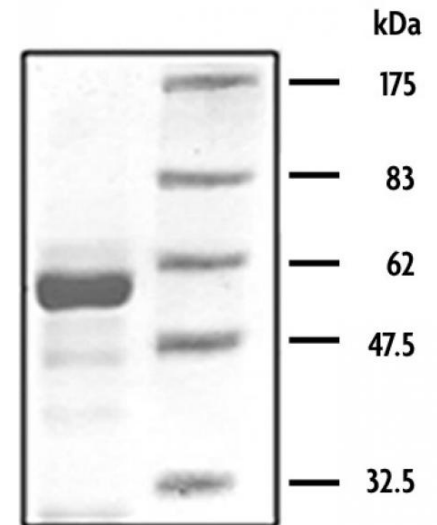
Background: SUV420H2 (Suppressor Of Variegation 4-20 Homolog 2 (Drosophila)) is a histone methyltransferase that specifically trimethylates Lys-20 of histone H4. H4 Lys -20 trimethylation represents a specific tag for epigenetic transcriptional repression. Mainly functions in pericentric heterochromatin regions, thereby playing a central role in the establishment of constitutive heterochromatin in these regions. SUV420H1 is targeted to histone H3 via its interaction with RB1 family proteins (RB1, RBL1 and RBL2) (By similarity).

Protein Details: SUV420H2 is a histone methyltransferase specific for trimethylation of Histone H4 at lysine 20. Amino acids 2-280 of SUV420H2 (accession number NM_032701) were expressed in *E. coli* with an N-terminal GST tag (MW= 58 kDa).

Application Notes: Recombinant SUV420H2 is suitable for use in histone methyltransferase (HMT) assays. It is useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



SUV420H2 protein gel.
SUV420H2 run on an SDS-PAGE gel and stained with Coomassie blue.