

Recombinant SUV420H1 protein

Catalog No: 31360

Expressed In: *E. coli*

Quantity: 50 µg

Concentration: 0.3 µg/µl

Source: Human

Buffer Contents: 50 µg recombinant SUV420H1 supplied at a concentration of 0.3 µg/µl in a buffer of 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, and 3 mM DTT.

Protein Details: SUV420H1 is a histone methyltransferase specific for dimethylation of Histone H4 at lysine 20. Amino acids 2-387 of human SUV420H1 (accession number NM_017635) were expressed in *E. coli* with an N-terminal GST tag (MW= 70 kDa).

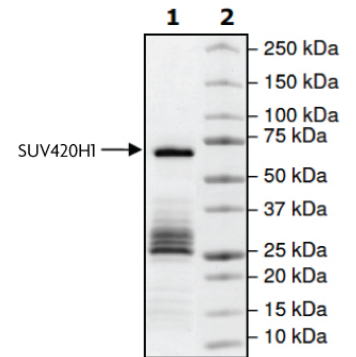
Application Notes: Recombinant SUV420H1 is suitable for use in histone methyltransferase (HMT) assays. This protein is useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Specific Activity: 0.0009 pmol/min/µg.

Assay conditions: Prepare a 50 µl reaction mix (50 mM TRIS pH 8.8, 5 mM MgCl₂, 4 mM DTT, 0.05% Tween-20, 800 µM S-adenosylmethionine, and SUV420H1) and add to substrate-coated wells. Incubate for 4 hr. Add primary antibody, incubate 1 hr. Next, incubate 30 minutes with the HRP-labeled secondary antibody followed by chemiluminescent detection.

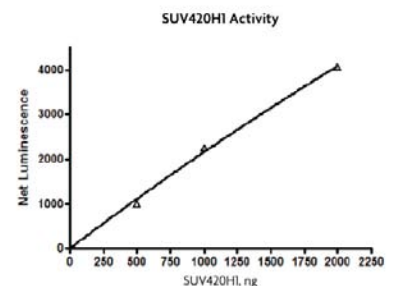
Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

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



SUV420H1 protein gel.

SUV420H1 run on a 4-20% SDS-PAGE gel and stained with Coomassie blue. Arrow indicates recombinant SUV420H1.



SUV420H1 activity assay.

Recombinant SUV420H1 activity measured using a fluorescent histone methyltransferase assay.