

Recombinant SETD3 protein

Catalog No: 81279, 81979

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.5 µg/µl

Source: Human

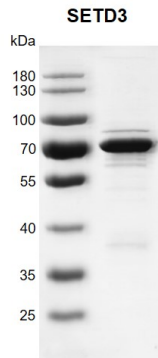
Buffer Contents: Recombinant SETD3 protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, and 0.5 mM TCEP.

Background: SETD3 (SET Domain Containing 3, Actin Histidine Methyltransferase), is a protein-histidine N-methyltransferase that specifically mediates methylation of actin at 'His-73'. Histidine methylation of actin is required for smooth muscle contraction of the laboring uterus during delivery. According to researches, SETD3 does not have proteinlysine N-methyltransferase activity and probably only catalyzes histidine methylation of actin.

Protein Details: Full length human SETD3 protein (accession number NP_115609.2 was expressed in a baculovirus expression system and contains an N-terminal FLAG-Tag. The molecular weight of SETD3 is 68.5 kDa.

Application Notes: Recombinant SETD3 protein is suitable for use in the study of enzyme kinetics, inhibitor screening, 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 arrival.

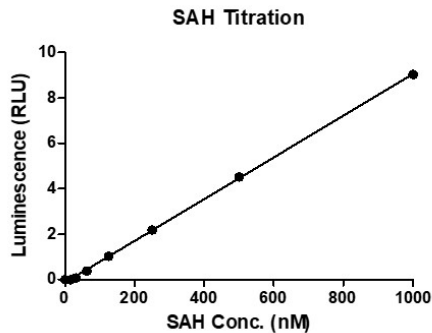


Recombinant SETD3 protein gel

10% SDS-PAGE Coomassie staining

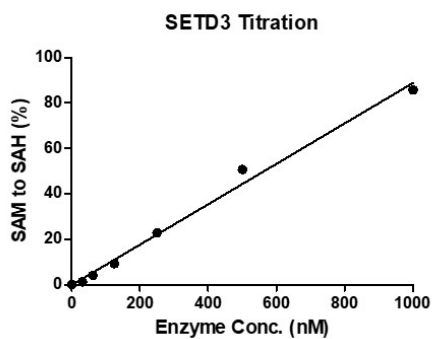
MW: 68.5 kDa

Purity: >90%



MTase-Glo Assay for Recombinant SETD3 protein methyltransferase activity

1 μ M beta-actin peptide (66-TLKYPIDHGVITNWD-80) and 1 μ M SAM were incubated with different concentrations of recombinant SETD3 protein in an 8 μ l reaction system containing 50 mM Tris-HCl pH 8.6, 0.02% Triton X-100, 2 mM MgCl₂, 1 mM TCEP for 1 hr. 5 \times MTase-Glo Reagent was added to the products and incubated for 30 min. Then MTase-Glo Detection was added and luminescence were read after another 30 min incubation. SAH standard curve (0-1 μ M) was performed following the same protocol.



MTase-Glo Assay for Recombinant SETD3 protein methyltransferase activity

1 μ M beta-actin peptide (66-TLKYPIDHGVITNWD-80) and 1 μ M SAM were incubated with different concentrations of recombinant SETD3 protein in an 8 μ l reaction system containing 50 mM Tris-HCl pH 8.6, 0.02% Triton X-100, 2 mM MgCl₂, 1 mM TCEP for 1 hr. 5 \times MTase-Glo Reagent was added to the products and incubated for 30 min. Then MTase-Glo Detection was added and luminescence were read after another 30 min incubation. SAH standard curve (0-1 μ M) was performed following the same protocol.