

Recombinant SETD6 protein

Catalog No: 81282, 81982

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.25 µg/µl

Source: Human

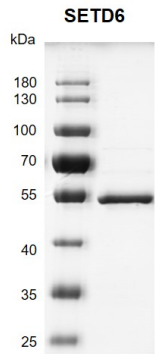
Buffer Contents: Recombinant SETD6 protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, and 0.5 mM TCEP.

Background: SETD6 (SET Domain Containing 6, Protein Lysine Methyltransferase) is a member of SETD family, which has methyltransferase activity. This protein was reported to monomethylate 'Lys-8' of histone H2AZ, which is involved in nuclear receptor-dependent transcription. SETD6 was also reported to monomethylates 'Lys-310' of the RELA subunit of NF kappa-B complex, leading to down-regulate NF-kappa-B transcription factor activity. It may interact with several endogenous proteins which are involved in nuclear hormone receptor signaling. However, the real substrates remain unknown for a very long time.

Protein Details: Full length human SETD6 protein (accession number NP_079136.2) was expressed in a baculovirus expression system and contains an N-terminal FLAG tag. The molecular weight of SETD6 is 52.1 kDa.

Application Notes: Recombinant SETD6 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.

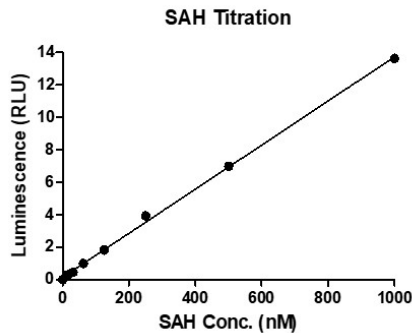


Recombinant SETD6 protein gel

10% SDS-PAGE with Coomassie staining.

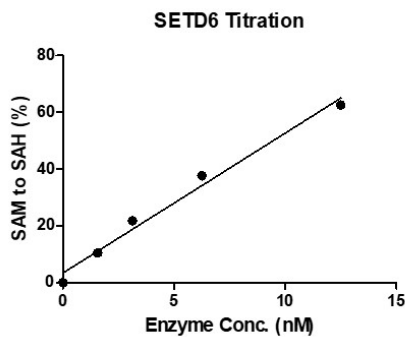
MW: 52.1 kDa

Purity: >90%



MTase-Glo Assay for Recombinant SETD6 protein methyltransferase activity

1 μ M substrate peptide and 1 μ M SAM were incubated with different concentrations of recombinant SETD6 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 was read after another 30 min incubation. SAH standard curve (0-1 μ M) was performed following the same protocol



MTase-Glo Assay for Recombinant SETD6 protein methyltransferase activity

1 μ M substrate peptide and 1 μ M SAM were incubated with different concentrations of recombinant SETD6 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 was read after another 30 min incubation. SAH standard curve (0-1 μ M) was performed following the same protocol