

Recombinant SETD4 protein

Catalog No: 81280, 81980

Expressed In: Baculovirus

Quantity: 20, 1000 µg

Concentration: 0.1 µg/µl

Source: Human

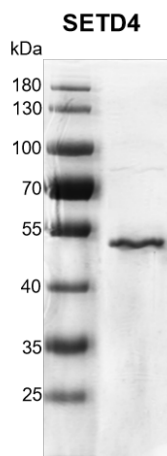
Buffer Contents: Recombinant SETD4 protein is supplied in 25 mM HEPES-NaOH pH 7.5, 300 mM NaCl, 10% glycerol, 0.04% Triton X-100, 0.5 mM TCEP.

Background: **SETD4** (SET domain-containing protein 4), Histone-lysine N-methyltransferase that acts as a regulator of cell proliferation, cell differentiation and inflammatory response. Regulates the inflammatory response by mediating mono and dimethylation of 'Lys-4' of histone H3 (H3K4me1 and H3K4me2, respectively), leading to activate the transcription of proinflammatory cytokines IL6 and TNF-alpha. Also involved in the regulation of stem cell quiescence by catalyzing the trimethylation of 'Lys-20' of histone H4 (H4K20me3), thereby promoting heterochromatin formation. Involved in proliferation, migration, paracrine and myogenic differentiation of bone marrow mesenchymal stem cells (BMSCs).

Protein Details: Full length human SETD4 protein (accession number NP_059134.1) was expressed in baculovirus expression system and contains an N-terminal FLAG tag. The molecular weight of SETD4 is 51.7 kDa.

Application Notes: Recombinant SETD4 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 for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.



Recombinant SETD4 protein

10% SDS-PAGE with Coomassie blue staining

MW: 51.7 kDa

Purity: >90%