

Recombinant METTL21D (VCPKMT) protein

Catalog No: 81210

Expressed In: Baculovirus

Quantity: 20 µg

Concentration: 0.15 µg/µl

Source: Human

Buffer Contents: Recombinant METTL21D (VCPKMT) 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.

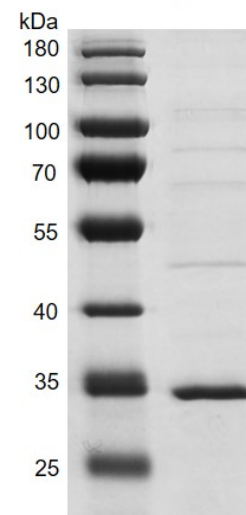
Background: METTL21D (Methyltransferase-Like Protein 21D), also known as VCPKMT (Valosin Containing Protein Lysine Methyltransferase), is a member of the Family 16 group of methyltransferases. It is responsible for the trimethylation of VCP (also called p97) on Lys 315. VCP is a highly conserved member of the AAA+-ATPases (ATPases associated with various activities) family, involved in several cellular pathways including protein degradation, endoplasmic reticulum-associated degradation, cell cycle regulation, membrane fusion and DNA damage response. Recently, it has been shown that VCPKMT is involved in cancer metastasis formation, but the physiological role of K315me3 in VCP has not been established and also the effect of K315me3 on the ATPase activity of VCP is not yet clear.

Protein Details: Recombinant METTL21D (VCPKMT) protein was expressed in a baculovirus expression system as the full length human protein (accession number NP_078834.2) with an N-terminal FLAG-Tag. The molecular weight of the protein is 27 kDa.

Application Notes: This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data this product is shown.

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.

METTL21D (VCPKMT)



Recombinant METTL21D (VCPKMT) protein gel

10% SDS-PAGE gel with Coomassie blue staining
MW: 27 kDa
Purity: >90%