

SETMAR Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # S347-30G

Lot # N343-3

Product Description

Recombinant full-length human SETMAR was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC011635](#).

Gene Aliases

METNASE

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

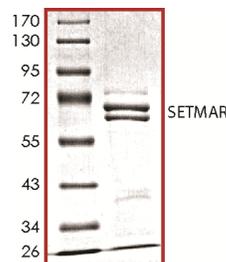
Scientific Background

SETMAR or SET domain and mariner transposase fusion gene is a nonhomologous end-joining repair protein that regulates genomic integration of exogenous DNA by opening chromatin and facilitating joining of DNA ends (1). SETMAR has histone methyltransferase activity and methylates 'Lys-4' and 'Lys-36' of histone H3. SETMAR also has DNA nicking activity and may play a role in DNA repair (2). Human Pso4 forms a stable complex with SETMAR that regulates Metnase function in DNA repair. SETMAR has sequence-specific DNA-binding activity and recognizes the 19-mer core of the 5'-terminal inverted repeats (TIRs) of the Hsmar1 element.

References

- Lee, S.-H.et.al: The SET domain protein Metnase mediates foreign DNA integration and links integration to nonhomologous end-joining repair. Proc. Nat. Acad. Sci. 102: 18075-18080, 2005.
- Liu, D.et.al: The human SETMAR protein preserves most of the activities of the ancestral Hsmar1 transposase. Molec. Cell. Biol. 27: 1125-1132, 2007.

Purity



The purity of SETMAR was determined to be **>85%** by densitometry.
Approx. MW **66-69 kDa**.

SETMAR Protein

Full length recombinant protein expressed in Sf9 cells

Catalog Number	S347-30G
Specific Lot Number	N343-3
Purity	>85%
Concentration	0.2 µg/µl
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.