

KAT8 (MYST1) Protein

Recombinant human protein expressed in Sf9 cells

Catalog # K317-31G

Lot # 1196-1

Product Description

Recombinant human KAT8 (2-467; contains the catalytic domain) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [BC037773](#).

Gene Aliases

KAT8, MYST1, hMOF, MOF, FLJ14040

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM 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.

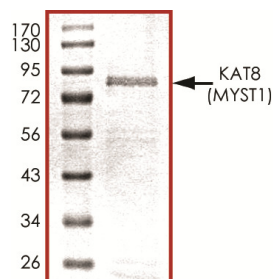
Scientific Background

KAT8 (MYST1) or MYST histone acetyltransferase 1 is a signaling protein that belongs to the MYST family of histone acetyl transferases (HATs) and was originally isolated as an HIV-1 TAT-interactive protein which play important roles in regulating chromatin remodeling, transcription and other nuclear processes by acetylating histone and nonhistone proteins. MYST1 have a chromodomain that involved in protein-protein interactions and targeting transcriptional regulators to chromatin (1). MYST histone acetyltransferase 1 is also involved in ATM function (2).

References

1. Neal, K. C. et.al: A new human member of the MYST family of histone acetyl transferases with high sequence similarity to Drosophila MOF. Biochim. Biophys. Acta 1490: 170-174, 2000.
2. Gupta, A. et.al: Involvement of human MOF in ATM function. Molec. Cell. Biol. 25: 5292-5305, 2005.

Purity



The purity was determined to be **>80%** by densitometry. Approx. MW **83 kDa**.

KAT8 (MYST1) Protein

Recombinant protein expressed in Sf9 cells

Catalog Number	K317-31G
Specific Lot Number	1196-1
Purity	>80%
Concentration	0.05 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.