

PRODUCT INFORMATION



TIP60 (human, recombinant)

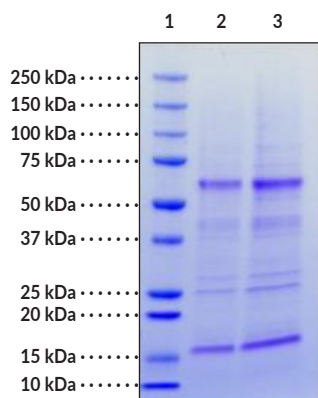
Item No. 10783

Overview and Properties

Synonyms:	cPLA(2) Interacting Protein, Esa1, Histone acetyltransferase KAT5, HIV-1 Tat Interacting Protein 60 kDa, Hs.6364, HTATIP, KAT5, PLIP
Source:	Recombinant N-terminal hexahistidine-tagged protein expressed in Sf21 cells
Amino acids:	2-513 (Full length, Isoform 2)
Uniprot No.:	Q92993
Molecular Weight:	60.3 kDa
Storage:	-80°C (as supplied); avoid freeze/thaw cycles by aliquoting protein
Stability:	≥6 months
Purity:	≥50% (estimated by SDS-PAGE)
Supplied in:	50 mM Tris-HCl, pH 7.5, with 300 mM sodium chloride and 10% glycerol
Protein Concentration:	<i>batch specific</i> mg/ml
Activity:	<i>batch specific</i> U/ml
Specific Activity:	<i>batch specific</i> U/mg
Unit Definition:	One unit is defined as the amount of enzyme required to produce 1 pmole of methylated peptide per minute at 37°C in 100 mM HEPES, pH 7.5, 0.8% Triton X-100 containing 100 μM histone H4 peptide (amino acids 1-21, Item No. 10854), and acetyl-CoA. Activity was determined using Cayman's HAT Inhibitor Screening Assay Kit (Item No. 10006515).

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

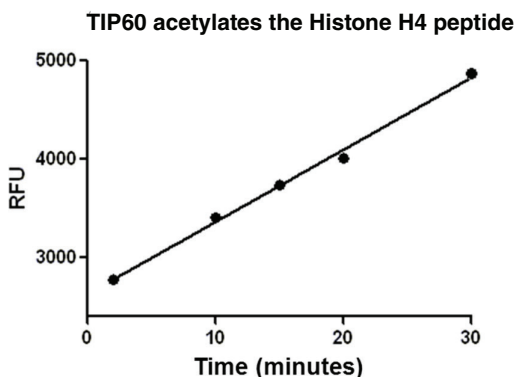
Images



Lane 1: MW Markers
Lane 2: TIP60 (2 μg)
Lane 3: TIP60 (4 μg)

SDS-PAGE Analysis of TIP60.

Representative gel image shown; actual purity may vary between each batch.



WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 08/24/2021

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

TIP60 is a member of the MYST family of lysine acetyl transferases. It has been shown to acetylate histones, p53, and the Ataxia Telangiectasia Mutant (ATM) protein kinase.¹ As a component of the TIP60-p400 complex, TIP60 acetylates transcription repressors in order to inhibit gene transcription.² In complexes without p400, TIP60 is part of a chromatin remodeling complex which plays an important role in double-stranded DNA break repair. Acetylation of histone γ H2AX allows formation of the γ H2AX repair foci, and acetylation of ATM is one of the early steps in double-stranded DNA break repair. Furthermore, TIP60 is bound by the cytoplasmic tail of β -amyloid precursor protein which drives gene transcription, apoptosis, and potentially a role in Alzheimer's disease pathogenesis.³

References

1. Sun, Y., Jiang, X., and Price, B.D. Tip60: Connecting chromatin to DNA damage signaling. *Cell Cycle* **9(5)**, 930-936 (2010).
2. Fazio, T.G., Hurr, J.T., and Panning, B. Chromatin regulation Tip(60)s the balance in embryonic stem cell self-renewal. *Cell Cycle* **7(21)**, 3302-3306 (2008).
3. Wu, J., Wang, J., Li, M., *et al.* Small molecule inhibitors of histone acetyltransferase Tip60. *Bioorg. Chem.* **39(1)**, 53-58 (2011).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM