

CERTIFICATE of ANALYSIS



TIP60 (human, recombinant)

Item No. 10783 • Batch No. 0805049 • Expires 21APR2026

Overview

Formulation:	50 mM Tris-HCl, pH 7.5, with 300 mM sodium chloride and 10% glycerol
Concentration:	1.0 mg/mL
Storage/stability:	-80°C/6mo
Specific activity:	5,038.00 U/mg
Unit definition:	One unit is defined as the amount of enzyme required to produce 1 pmole of acetylated peptide per minute at 37°C in 100 mM HEPES, pH 7.5, 0.8% Triton X-100 containing 100 μ M 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).
Activity:	5,038.00 U/ml
Purity:	$\geq$ 50%

Reviewed and approved by: Beth Hurst

Approved date: 22OCT2025

Elizabeth Hurst

WARNING

THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. IT IS THE RESPONSIBILITY OF THE PURCHASER TO DETERMINE SUITABILITY FOR OTHER APPLICATIONS.

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, 10/12/2018

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