

CERTIFICATE of ANALYSIS



HDAC8 (human, recombinant)

Item No. 19380 • Batch No. 0683705 • Expires 10JAN2026

Overview

Formulation:	50 mM Tris, pH 7.5, with 200 mM NaCl, 2.5 mM potassium chloride, 1 mM TCEP, and 25% glycerol
Storage/stability:	-80°C/6mo
Specific activity:	5.16 U/mg
Unit definition:	One unit is the amount of enzyme required to release 1 nmol/min acetate from 100 μ M acetylated p53 peptide (Item No. 10010995) at 37°C in 25 mM Tris, pH 8.0, 137 mM sodium chloride, 2.7 mM KCl, and 1 mM MgCl ₂ .
Activity:	2.17 U/ml
Purity:	≥80% estimated by SDS-PAGE

Reviewed and approved by: Carly Norris

Approved date: 02AUG2023

A handwritten signature in blue ink that reads "Carly Norris".

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