

CERTIFICATE of ANALYSIS



SIRT3 (human, recombinant)

Item No. 10011194 • Batch No. 0807039 • Expires 03JUN2027

Overview

Formulation:	50 mM sodium phosphate, pH 7.2, with 100 mM sodium chloride and 20% glycerol
Concentration:	0.32 mg/mL
Storage/stability:	-80°C/1yr
Specific activity:	27.80 U/mg
Unit definition:	One unit is defined as the amount of enzyme required to produce 1 nmole of 7-amino-4-methylcoumarin per minute at 25°C in 50 mM Tris-HCl, pH 8.0, containing 137 mM NaCl, 2.7 mM KCl, 1 mM MgCl ₂ , 125 µM p53 amino acids 317-320 (Gln-Pro-Lys-Lys(e-acetyl)-AMC), and 6 mM NAD ⁺
Activity:	8.90 U/ml
Purity:	≥60% estimated by SDS-PAGE

Reviewed and approved by: Beth Hurst

Approved date: 03JUN2026

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