

CERTIFICATE of ANALYSIS



Histone H3K4Me3 Monoclonal Antibody (RM340)

Item No. 32130 • Batch No. 0819232 • Expires 16MAR2027

Overview

Formulation:	100 µg of protein A-affinity purified monoclonal antibody
Concentration:	1 mg/mL
Storage/stability:	-20°C/1yr
Applications:	ChIP, ELISA, multiplex-based assays
Cross reactivity:	(-)H3K4Me1(-)H3K4Me2(-)H3K9Me1(-)H3K9Me2(-)H3K9Me3(-)H3K27Me1(-)H3K27Me2(-)H3K27Me3(-)H3K36Me1(-)H3K36Me2(-)H3K36Me3(-)Unmodified H3K4(-)H4K20Me1(-)H4K20Me2(-)H4K20Me3
Species reactivity:	(+)Vertebrates(+)H3K4Me3

Reviewed and approved by: Justine Albertson

Approved date: 13MAR2026

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