

PRODUCT INFORMATION

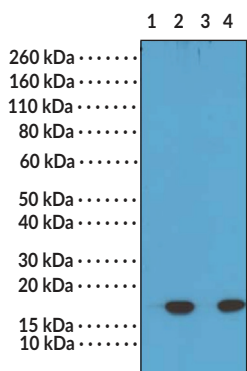
Histone H3^{K36M} Rabbit Monoclonal Antibody (Clone RM193)

Item No. 31547

Overview and Properties

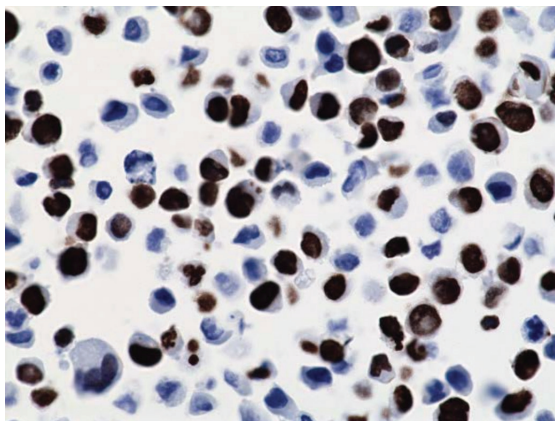
Contents:	This vial contains 100 or 400 µl of protein A-affinity purified monoclonal antibody.
Immunogen:	A peptide corresponding to histone H3K36M mutant
Cross Reactivity:	(-) Wild-type histone H3
Species Reactivity:	(+) Histone H3K36M mutant
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RM193
Host:	Rabbit
Isotype:	IgG
Applications:	ELISA, immunohistochemistry (IHC), and Western blot (WB); the recommended starting dilution for ELISA is 1:250-1:2,500, 1:100-1:500 for IHC, and 1:200-1:1,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Lane 1: H3.1 WT
Lane 2: H3.1 K36M
Lane 3: H3.3 WT
Lane 4: H3.3 K36M

WB analysis of cell lysates prepared from 293T transfected with a DNA construct encoding wild-type or K36M mutant proteins of histone H3.1 and H3.3, using Histone H3^{K36M} Rabbit Monoclonal Antibody (Clone RM193).



Immunohistochemical staining of formalin-fixed and paraffin-embedded 293T cells transfected with a DNA construct encoding histone H3K36M, stained with Histone H3^{K36M} Rabbit Monoclonal Antibody (Clone RM193).

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, 02/25/2025

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

Histone H3 undergoes many modifications including acetylation, methylation, and phosphorylation that are important for regulation of gene transcription. Lysine 36 of histone 3 (H3K36) has functionally relevant acetylation and methylation states with roles in transcriptional activation. More than 90% of chondroblastomas contain a heterozygous mutation, replacing H3K36 with methionine 36 (H3K36M) in the histone H3 variant H3.3.¹ This antibody reacts to the H3K36M mutant and does not detect wild-type histone H3.

Reference

1. Fang, D., Gan, H., Lee, J.-H., *et al.* The histone H3.3K36M mutation reprograms the epigenome of chondroblastomas. *Science* **352**(6291), 1344-1348 (2016).

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