

CERTIFICATE of ANALYSIS



Anti-Histone H3 pan Rabbit Monoclonal Antibody (Clone RM188)

Item No. 20724 • Batch No. 0700186 • Expires 18JAN2025

Overview

Formulation:	100 µg of protein A affinity-purified antibody from an animal origin-free culture supernatant 50% Glycerol/PBS with 1% BSA and 0.09% sodium azide
Concentration:	1.0 mg/mL
Storage/Stability:	-20°C/1yr
Applications:	ELISA, ICC, multiplex-based assays, WB
Cross reactivity:	(+)All histone H3 proteins, independent of post-translational modifications, (-)H2A, (-)H2B, (-)H4
Species reactivity:	(Independent)Species

Reviewed and approved by: Emily Tucker

Approved date: 19JAN2024

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