

CERTIFICATE of ANALYSIS



4-hydroxy Nonenal Monoclonal Antibody (HNEJ-1/Clone IG10) - HRP Conjugated
Item No. 45025 • Batch No. 0818915 • Expires 05MAR2027

Overview

Formulation:	100 µg protein G purified antibody in PBS with 50% glycerol, 0.05% ProClin300 and proprietary stabilizers
Concentration:	0.5 mg/mL
Storage/stability:	-20°C/2yr
Applications:	WB

Reviewed and approved by: Beth Hurst

Approved date: 06MAR2026

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