

CERTIFICATE of ANALYSIS



MR1 Monoclonal Antibody (Clone 2H7)

Item No. 34257 • Batch No. 0711255 • Expires 22MAY2027

Overview

Formulation: 5 mg of protein G-affinity purified monoclonal antibody in PBS
Concentration: 2.65 mg/mL
Storage/Stability: -20°C/3yr
Applications: ELISA, FC, IP (mass spec confirmed), WB
Cross reactivity: (+)MR1
Species reactivity: (+)Human

Reviewed and approved by: Jeremy Bickel

Approved date: 23MAY2024

A handwritten signature in black ink, appearing to read "Jeremy Bickel", written over a light blue horizontal line.

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