

CERTIFICATE of ANALYSIS



M13 Monoclonal Antibody (Clone MM05)

Item No. 38076 • Batch No. 0822951 • Expires 19JUN2027

Overview

Formulation:	50, 100, or 200 µl of protein A-affinity purified recombinant monoclonal antibody in 0.2 µm filtered solution in PBS with 0.02% Tween 80, pH 7.4
Concentration:	1 mg/mL *Buffer
Storage/stability:	-80°C/1yr
Applications:	ELISA
Cross reactivity:	(+)Phage coat proteins
Species reactivity:	(+)M13 bacteriophage(+)fd bacteriophage

Reviewed and approved by: Justine Albertson

Approved date: 10JUN2026



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