

CERTIFICATE of ANALYSIS



AcMNPV Major Envelope Glycoprotein Neutralizing Antibody (Clone M001)

Item No. 37014 • Batch No. 0713400 • Expires 17JUN2025

Overview

Formulation:	200 or 500 µg of protein A-affinity purified monoclonal antibody 0.2 µm filtered solution in Histidine and Arginine buffer containing 120mM NaCl, 0.02% Tween 80, pH6.0
Concentration:	10.02 mg/mL
Storage/Stability:	-80°C/1yr
Applications:	MN
Cross reactivity:	(+)Major envelope glycoprotein
Species reactivity:	(+)AcMNPV

Reviewed and approved by: Emily Tucker

Approved date: 13JUN2024

A handwritten signature in black ink, appearing to read "Emily Tucker", with a stylized flourish at the end.

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