

CERTIFICATE of ANALYSIS



SARS-CoV/SARS-CoV-2 Nucleocapsid Protein Monoclonal Antibody (Clone 08)

Item No. 31984 • Batch No. 0719579 • Expires 26AUG2025

Overview

Formulation:	50 or 100 µl of protein A-affinity purified monoclonal antibody in PBS
Concentration:	1.24 mg/mL
Storage/Stability:	-80°C/1yr
Applications:	ELISA, FC, IHC-P, WB
Cross reactivity:	(+)Nucleocapsid protein, (+)SARS-CoV-2 Alpha (B.1.1.7) nucleocapsid protein (D3L, R203K, G204R, S235F), (+)SARS-CoV-2 Alpha (B.1.1.7) nucleocapsid protein (D3L, S235F), (+)SARS-CoV-2 Alpha (B.1.1.7) nucleocapsid protein (I292T), (+)SARS-CoV-2 Alpha (B.1.1.7) nucleocapsid protein (R203K, G204R), (+)SARS-CoV-2 Alpha (B.1.1.7/B.1.237) nucleocapsid protein (S194L), (+)SARS-CoV-2 Alpha/Beta (B.1.1.7/B.1.351/A.2.2) nucleocapsid protein (P13L), (+)SARS-CoV-2 Beta (B.1.351/B.1.351.2/B.1.351.3/B.1.427/B.1.429) nucleocapsid protein (T205I), (+)SARS-CoV-2 Delta (B.1.617) nucleocapsid protein (D377Y), (+)SARS-CoV-2 Delta (B.1.617) nucleocapsid protein (R230M, D377Y), (+)SARS-CoV-2 Delta (B.1.617.3) nucleocapsid protein (P67S, R203M, D377Y), (+)SARS-CoV-2 Eta (B.1.525) nucleocapsid protein (A12G, T205I), (+)SARS-CoV-2 Gamma (P.1/P.1.1/P.1.2) nucleocapsid protein (P80R)
Species reactivity:	(+)SARS-CoV, (+)SARS-CoV-2, (+)SARS-CoV-2 Delta (B.1.617), (+)SARS-CoV-2 Delta (B.1.617.2), (+)SARS-CoV-2 Omicron (B.1.1.529), (+)SARS-CoV-2 Omicron (BA.2), (+)SARS-CoV-2 Omicron (BA.4), (+)SARS-CoV-2 Omicron XE (BA.1 x BA.2), (-)HCoV-229E, (-)HCoV-HKU1 (isolate N5), (-)HCoV-NL63, (-)HCoV-OC43, (-)MERS-CoV

Reviewed and approved by: Emily Tucker

Approved date: 26AUG2024

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