

CERTIFICATE of ANALYSIS



SARS-CoV/SARS-CoV-2 Spike Glycoprotein RBD Chimeric Monoclonal Antibody (Clone D005)

Item No. 30832 • Batch No. 0617636 • Expires 01JUN2022

Overview

Formulation: 50 µl of protein A-purified monoclonal antibody

Storage/Stability: -80°C/1yr

Applications: ELISA

Cross reactivity: (+)SARS-CoV spike glycoprotein S1 subunit, (+)SARS-CoV-2 (BA.2.75) spike glycoprotein S1+S2 trimer, (+)SARS-CoV-2 (BA.4.6) spike glycoprotein S1+S2 trimer, (+)SARS-CoV-2 Delta (B.1.617.2) spike glycoprotein S1 subunit, (+)SARS-CoV-2 Delta (B.1.617.2) spike glycoprotein S1+S2, (+)SARS-CoV-2 Delta (B.1.617.2) spike glycoprotein S1+S2 trimer, (+)SARS-CoV-2 Omicron (B.1.1.529) spike glycoprotein S1 subunit, (+)SARS-CoV-2 Omicron (B.1.1.529) spike glycoprotein S1+S2 trimer, (+)SARS-CoV-2 Omicron (BA.1.1) spike glycoprotein S1+S2 trimer, (+)SARS-CoV-2 Omicron (BA.2) spike glycoprotein S1 subunit, (+)SARS-CoV-2 Omicron (BA.2) spike glycoprotein S1+S2 trimer, (+)SARS-CoV-2 Omicron (BA.2.75.2) spike glycoprotein S1+S2 trimer, (+)SARS-CoV-2 Omicron (BF.7) spike glycoprotein S1+S2 trimer, (+)SARS-CoV-2 Omicron (BQ.1.1) spike glycoprotein S1+S2 trimer, (+)SARS-CoV-2 XD (BA.1 x AY.4) spike glycoprotein S1+S2 trimer, (+)SARS-CoV-2 spike glycoprotein S1 subunit, (-)HCoV-229E spike glycoprotein S1 subunit, (-)HCoV-HKU1 (isolate N1) spike glycoprotein S1 subunit, (-)HCoV-HKU1 (isolate N5) spike glycoprotein S1 subunit, (-)HCoV-NL63 spike glycoprotein S1 subunit, (-)HCoV-OC43 spike glycoprotein S1+S2 ECD, (-)MERS-CoV spike glycoprotein S1 subunit, (-)SARS-CoV-2 Delta (B.1.617.2) spike glycoprotein S1 subunit NTD, (-)SARS-CoV-2 Omicron (BA.2) spike glycoprotein S1 subunit NTD

Species reactivity: (+)SARS-CoV, (+)SARS-CoV-2, (+)SARS-CoV-2 (BA.2.3.20), (+)SARS-CoV-2 (BA.2.75), (+)SARS-CoV-2 Delta (B.1.617.2), (+)SARS-CoV-2 Omicron (B.1.1.529), (+)SARS-CoV-2 Omicron (BA.1.1), (+)SARS-CoV-2 Omicron (BA.2), (+)SARS-CoV-2 Omicron (BA.2.12.1), (+)SARS-CoV-2 Omicron (BA.2.75.2), (+)SARS-CoV-2 Omicron (BA.4), (+)SARS-CoV-2 Omicron (BA.4.6/BF.7), (+)SARS-CoV-2 Omicron (BA.5), (+)SARS-CoV-2 Omicron (BQ.1.1), (+)SARS-CoV-2 Omicron (XBB)

Reviewed and approved by: Kyle Bale

Approved date: 01JUN2021

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