

CERTIFICATE of ANALYSIS



ACE2 (human) Monoclonal Antibody (Clone AC384; preservative free)

Item No. 32028 • Batch No. 0680366 • Expires 20JUN2024

Overview

Formulation:	100 or 500 µg of Protein G-purified monoclonal antibody, stored in PBS.
Concentration:	1 mg/ml
Storage/Stability:	-20°C/1yr
Applications:	ELISA (direct or indirect), WB (using ECL)
Species reactivity:	(+)Human
Biofunctional assay:	Blocks the binding of human ACE2 to the Spike protein of SARS-Cov-2, but does not affect ACE2 activity. Optimal conditions should be determined individually for each application.

Reviewed and approved by: Judy Hines

Approved date: 21JUN2023

A handwritten signature in black ink, appearing to read "Judy P. Hines", 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