



Extension Memo

Page 1 of 1

Product Description: ACE2 Calibration Reagent
Catalog Number(s): C01ADG-2
Lot Number: A0080309
New Expiration Date: 30 Apr 2026

Based on additional testing, the original date of expiration for this calibrator has been extended by 12 months.

Function	Name	Signature	Date
Quality	Karen Hamilla		10 Mar 2025



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: ACE2 Calibration Reagent

Catalog Number: C01ADG-2

Lot Number: A0080309

Expiration Date: 30 APR 2025


Storage Temperature: 2-8°C

Concentration Assignment:

	Stock Concentration	Unit of Measure
ACE2 Calibration Reagent	167	Units/mL

Dilution to Top of Curve: 10X

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Carissa B. Hubbard	Quality		23 DEC 2021