



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: ACE2 Calibration Reagent

Catalog Number: C01ADG-2

Lot Number: A0080301

Expiration Date: 31 MAR 2025

Storage Temperature: 2-8°C

Concentration Assignment:

	Stock Concentration	Unit of Measure
ACE2 Calibration Reagent	192	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	Karen Hamilla	Quality	<i>Karen Hamilla</i>	06 OCT 2021