



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: ACE2 Calibration Reagent 2

Catalog Number: C01AJZ-2

Lot Number: A0080327

Expiration Date: 31 OCT 2025


Storage Temperature: 2-8°C

Concentration Assignment:

	Stock Concentration	Unit of Measure
ACE2 Calibration Reagent 2	217	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		10 MAY 2022