



CERTIFICATE OF ANALYSIS

Product Description: ACE2 Calibration Reagent
Catalog Number: C01ADG-2
Lot Number: A0080457
Expiration Date: 30 SEP 2029
Storage Temperature: 2-8°C

Concentration Assignment

	Stock Concentration	Unit of Measure
ACE2 Calibration Reagent	195	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		07 MAY 2026