

**CERTIFICATE OF ANALYSIS**


Product Description: ACE2 Calibration Reagent 2
Catalog Number: C01AJZ-2
Lot Number: A0080446
Expiration Date: 30 JUN 2029
Storage Temperature: 2-8°C

Concentration Assignment

	Stock Concentration	Unit of Measure
ACE2 Calibration Reagent 2	254	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		22 JAN 2026