



CERTIFICATE OF ANALYSIS

Product Description: Rat Clusterin Calibrator
Catalog Number: C03MG-2
Lot Number: A0040289
Expiration Date: 31 DEC 2026
Storage Temperature: ≤ -70°C

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
Clusterin	2220	ng/mL

Dilution to Top of Curve: 20 X

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		15 AUG 2024