



CERTIFICATE OF ANALYSIS

Product Description: Rat Clusterin Kit
Lot Number: K0040562
Expiration Date: 31 JUL 2026
Catalog Number: K153MGC-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-rat Clusterin Antibody	D0041042	2-8°C	31 JUL 2027
Rat Clusterin Plate	Z0057437	2-8°C	31 JUL 2026
Diluent 29	M0290036	≤ -10°C	30 SEP 2026
Rat Clusterin Calibrator	A0040285	≤ -70°C	30 SEP 2026

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

Spot	Name	Calibrator		Rat Clusterin Control 1		Rat Clusterin Control 2		Rat Clusterin Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Clusterin	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	Clusterin	2060	ng/mL

Dilution to top of Curve: 20X

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		30 MAY 2024