

**CERTIFICATE OF ANALYSIS**

**Product Description:** Rat Clusterin Kit  
**Lot Number:** K0040574  
**Expiration Date:** 30 SEP 2026  
**Catalog Number:** K153MGC-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-rat Clusterin Antibody	D0041062	2-8°C	31 DEC 2027
Rat Clusterin Plate	Z0057484	2-8°C	31 DEC 2026
Diluent 29	M0290036	≤ -10°C	30 SEP 2026
Rat Clusterin Calibrator	A0040289	≤ -70°C	31 DEC 2026

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

**Precision, Accuracy, and Sensitivity Specifications:**


		Calibrator		Rat Clusterin Control 1		Rat Clusterin Control 2		Rat Clusterin Control 3	
Spot	Name	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	2220	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
<b>Review/Approval</b>	Karen Hamilla	Quality		08 MAY 2025