

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Rat Clusterin Kit
Kit Lot Number: K0040563
Expiration Date: 31 Jan 2025
Kit Catalog Numbers: K153MGC-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Rat Clusterin Plate	Z0057421	2-8°C	31 May 2026
Anti-rat ClusterinAntibody	D0040977	2-8°C	30 Nov 2025
Rat Clusterin Calibrator	A0040264	≤ -70°C	31 Jan 2025
Diluent 29	M0290033	≤ -10°C	30 Nov 2025

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

Precision and Accuracy 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	2080	ng/mL

Dilution to Top of Curve = 20X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		04 Apr 2024