



**RFE MEMO**

Page 1 of 1

**Product Description(s):** Rat Clusterin Kit  
**Catalog Number(s) / Size(s):** K153MGC - Series  
**Lot Number:** K0040504

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0040504 has been extended to 31 MAY 2022.

The expiration date for the following kit component has been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Rat Clusterin Calibrator	A0040221	31 MAY 2021	31 MAY 2022

Function	Name	Signature	Date
Quality	Karen Hamilla		24 MAR 2021

This memo supersedes all other statements regarding date of expiration for this product.

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Rat Clusterin Kit  
**Kit Lot Number:** K0040504  
**Expiration Date:** 31 May 2021  
**Kit Catalog Numbers:** K153MGC-Series

**Precision and Accuracy 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	1998	ng/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Rat Clusterin Plate	Z0056770	2-8°C	31 Jan 2023
Anti-rat Clusterin Antibody	D0040934	2-8°C	29 Feb 2024
Rat Clusterin Calibrator	A0040221	≤ -70°C	31 May 2021
Diluent 29	M0290027	≤ -10°C	31 Dec 2022

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		05 FEB 2021