



**RFE MEMO**


Page 1 of 1

**Product Description(s):** Rat Clusterin Kit  
**Catalog Number(s) / Size(s):** K153MGC-Series  
**Lot Number:** K0040448

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0040448 has been extended to 31 OCT 2020.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Diluent 29	M0290023	31 OCT 2019	31 OCT 2020

Function	Name	Signature	Date
Quality	Karen Hamilla		14 AUG 2019

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

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** Rat Clusterin Kit  
**Kit Catalog Numbers:** K153MGC - Series  
**Kit Lot Number:** K0040448  
**Expiration Date:** 31 Oct 2019

### Acceptance Data/ Percent Recovery 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 Material Tested :

Spot	Name	Item	Conc	Units
1	Clusterin	20X Calibrator	1998	ng/mL
		Rat Clusterin Control 1	123	ng/mL
		Rat Clusterin Control 2	29.3	ng/mL
		Rat Clusterin Control 3	6.93	ng/mL

Concentrations provided in this table refer to the stock concentration of the calibrators. For top of the standard curve concentrations, please refer to the Standard Curve Data

### Standard Curve Data:

Clusterin - Spot 1		
Conc. ng/mL	ECL Signal	%CV
99.9	296,000	2.2
25.0	155,000	3.8
6.24	44,900	2.8
1.56	10,500	3.5
0.390	2,470	4.8
0.0976	694	2.7
0.0244	259	2.0
0.00	111	3.8

### CERTIFICATE OF ANALYSIS

Page 2 of 2

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Rat Clusterin Plate	Z0056192	2-8°C	28 Feb 2021
Anti-rat Clusterin Antibody	D0040802	2-8°C	28 Feb 2022
Rat Clusterin Calibrator	A0040221	≤ -70°C	31 May 2021
Blocker A Solution	Not Kit Specific	Not Kit Specific	N/A
Diluent 29	M0290023	≤ -10°C	31 Oct 2019
Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

**Statement:**

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	28 FEB 2019