



RFE MEMO

Page 1 of 1

Product Description(s): Rat Clusterin Kit
Catalog Number(s) / Size(s): K153MGC - Series
Lot Number: K0040399

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0040399 has been extended to 31 Mar 2019.

The expiration dates for the following kit components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Rat Clusterin Calibrator	A0040143	31 Mar 2018	31 Mar 2019
Diluent 29	M0290021	31 May 2018	31 Mar 2019

Function	Name	Signature	Date
Quality Assurance	Dawn Stollar		09 APR 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: K0040399
Expiration Date: 31 Mar 2018

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	Rat Clusterin	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Rat Clusterin	20X Calibrator	2160	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

Control data was generated utilizing internal reference controls.

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:

Rat Clusterin - Spot 1		
Conc. ng/mL	ECL Signal	%CV
108	215,000	2.3
27.0	137,000	3.4
6.75	47,300	2.2
1.69	11,900	2.4
0.422	2,860	6.0
0.105	751	3.3
0.0264	233	3.0
0.00	71	5.8

CERTIFICATE OF ANALYSIS

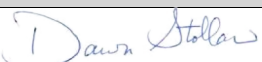
Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Rat Clusterin Plate	Z0055764	2-8°C	31 May 2019
Anti-rat Clusterin Anitibody	D0040657	2-8°C	30 Apr 2019
Rat Clusterin Calibrator	A0040143	≤ -70°C	31 Mar 2018
Blocker A Solution	Not Kit Specific	Not Kit Specific	N/A
Diluent 29	M0290021	≤ -10°C	31 May 2018
Diluent 29	M0290021	≤ -10°C	31 May 2018
MSD 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
Review/Approval	Dawn Stollar	Quality Assurance		09 APR 2019