

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Rat Clusterin Kit
Kit Catalog Numbers: K153MGC-Series
Kit Lot Number: K0040444
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	rt Clusterin	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	rt Clusterin	20X Calibrator	2080	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:

rt Clusterin - Spot 1		
Conc. ng/mL	ECL Signal	%CV
104	304,000	1.8
26.0	167,000	8.1
6.50	51,800	3.7
1.63	11,800	2.7
0.406	2,870	2.3
0.102	772	3.3
0.0254	274	2.3
0.00	104	3.1

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	A0040209	≤ -70°C	31 Aug 2020
Blocker A	Not Kit Specific	Not Kit Specific	N/A
Diluent 29	M0290023	≤ -10°C	31 Oct 2019
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		25 SEP 2018