



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Rat Clusterin Calibrator

**Catalog Number:** C03MG-2

**Lot Number:** A0040221

**Expiration Date:** 31 MAY 2021

**Storage Temperature:**  $\leq -70^{\circ}\text{C}$

### Concentration of Materials Tested:

	Assay
	Clusterin
Calibrator (ng/mL)	1998

**Concentration:** 20X

**Statement:** This 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>	16 JAN 2019