



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Rat B2M Calibrator
Catalog Number: C03KA-2
Lot Number: A0040192
Expiration Date: 31 AUG 2019
Storage Temperature: ≤ -70°C

Concentration of Materials Tested:

	Assay
	Rat B2M
Calibrator (ng/mL)	524

Statement: This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	28 APR 2017