



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Rat B2M Calibrator

Catalog Number: C03KA-2

Lot Number: A0040237

Expiration Date: 31 JAN 2022


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

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
Rat B2M	508	ng/mL

Dilution to Top of Curve: 20X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		09 JUN 2021