



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Rat B2M Calibrator

Catalog Number: C03KA-2

Lot Number: A0040275

Expiration Date: 30 APR 2025


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

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
B2M	528	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		03 JAN 2023