



CERTIFICATE OF ANALYSIS

Product Description: Rat B2M Calibrator
Catalog Number: C03KA-2
Lot Number: A0040303
Expiration Date: 31 OCT 2028
Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
B2M	578	ng/mL

Dilution to Top of Curve: 20 X

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		18 JUN 2026