



CERTIFICATE OF ANALYSIS

Product Description: Rat B2M Kit
Lot Number: K0040578
Expiration Date: 30 APR 2025
Catalog Number: K153KAC-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-rat B2M Antibody	D0041011	2-8°C	30 JUN 2026
Anti-rat B2M Biotinylated Capture Antibody	D0041010	2-8°C	30 JUN 2026
MSD GOLD™ 96-well Streptavidin Plate	Z0022275	2-8°C	30 SEP 2026
Diluent 5	M0050073	≤ -10°C	31 DEC 2026
Rat B2M Calibrator	A0040275	≤ -70°C	30 APR 2025

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

		Calibrator		Rat B2M Control 1		Rat B2M Control 2		Rat B2M Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	B2M	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%


		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	B2M	0.0488	80% - 120%	< 20%	20	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	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		15 OCT 2024