

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Rat B2M Kit
Kit Lot Number: K0040543
Expiration Date: 31 Jan 2025
Kit Catalog Numbers: K153KAC-Series

Precision and Accuracy 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%	20.0	80% - 120%	< 20%

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	B2M	528	ng/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
MSD GOLD™96-Well Streptavidin Plate	Z0021960	2-8°C	28 Feb 2025
Anti-rat B2M Antibody	D0041011	2-8°C	30 Jun 2026
Rat B2M Calibrator	A0040275	≤ -70°C	30 Apr 2025
Diluent 5	M0050071	≤ -10°C	31 Jan 2025
Anti-rat B2M Biotinylated Capture Antibody	D0041010	2-8°C	30 Jun 2026

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		30 JAN 2023