



Meso Scale Diagnostics, LLC.

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Rat B2M Kit
Kit Lot Number: K0040517
Expiration Date: 31 Jan 2022
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	508	ng/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
MSD GOLD 96-Well Streptavidin Plate	Z0021769	2-8°C	31 Aug 2023
Anti-rat B2M Antibody	D0040953	2-8°C	30 Nov 2024
Rat B2M Calibrator	A0040237	≤ -70°C	31 Jan 2022
Diluent 5	M0050068	≤ -10°C	31 Jan 2022
Anti-rat B2M Biotinylated Capture Antibody	D0040952	2-8°C	31 Dec 2024

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		09 AUG 2021