

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Rat B2M Kit
Kit Catalog Numbers: K153KAC-Series
Kit Lot Number: K0040471
Expiration Date: 30 Nov 2020

Acceptance Data/ Percent Recovery Specifications:

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

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	B2M	0.0488	80% - 120%	< 20%	20.0	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	B2M	20X Calibrator	516	ng/mL
		Rat B2M Control 1	4.79	ng/mL
		Rat B2M Control 2	1.1	ng/mL
		Rat B2M Control 3	0.279	ng/mL

Controls tested may not be the same lot as shipped.

Concentrations provided in this table refer to the stock concentration of the calibrators. For top of the standard curve concentrations, please refer to the Standard Curve Data

Standard Curve Data:

B2M - Spot 1		
Conc. ng/mL	ECL Signal	%CV
25.8	365,000	2.3
6.45	87,000	3.7
1.61	18,100	3.3
0.403	4,420	3.0
0.101	1,420	2.4
0.0252	652	2.8
0.00630	429	5.1
0.00	401	2.9

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
MSD GOLD™ 96-well Streptavidin Plate	Z0021434	2-8°C	30 Nov 2020
Anti-rat B2M Antibody	D0040850	2-8°C	30 Nov 2022
Anti-rat B2M Biotinylated Capture Antibody	D0040849	2-8°C	31 Dec 2022
Rat B2M Calibrator	A0040236	≤ -70°C	31 Oct 2021
Diluent 5	M0050067	≤ -10°C	30 Nov 2021
Blocker A Kit	Not Kit Specific	Not Kit Specific	N/A
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Rat B2M Control 1	Not Kit Specific
Rat B2M Control 2	Not Kit Specific
Rat B2M Control 3	Not Kit Specific

Statement:

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		08 AUG 2019