

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human MIP-1 β Kit

Kit Catalog Numbers: K151NR-Series, K156NR-Series

Kit Lot Number: K0081426

Expiration Date: 31 Jan 2022

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Chemokine Panel 1 (human) Control 1		Chemokine Panel 1 (human) Control 2		Chemokine Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	MIP-1 β	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	MIP-1 β	1.02	80% - 120%	< 20%	750	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	MIP-1 β	Calibrator	1330	pg/mL
		Chemokine Panel 1 (human) Control 1	1030	pg/mL
		Chemokine Panel 1 (human) Control 2	176	pg/mL
		Chemokine Panel 1 (human) Control 3	27.8	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

MIP-1 β - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1330	379,000	5.2
333	69,600	5.4
83.1	11,400	1.9
20.8	2,030	3.0
5.20	502	2.5
1.30	167	4.3
0.325	89	6.7
0.00	66	4.6

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 β Plate	Z0047262	2-8°C	30 Jun 2022
Anti-hu MIP-1 β Antibody	D0081121	2-8°C	31 Jan 2022
Chemokine Panel 1 (human) Calibrator Blend	A0080228	2-8°C	31 Aug 2024
Diluent 43	M0430036	\leq -10°C	31 May 2022
Diluent 3	M0030094	\leq -10°C	31 May 2022
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Plate Seals	Not Kit Specific
Wash Buffer (20X), 100mL	Not Kit Specific
Chemokine Panel 1 (human) Control 1	Not Kit Specific
Chemokine Panel 1 (human) Control 2	Not Kit Specific
Chemokine Panel 1 (human) 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		18 MAR 2020