

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: K0081471
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	356,000	4.9
333	58,200	5.9
83.1	8,170	2.5
20.8	1,120	2.6
5.20	241	2.5
1.30	102	7.0
0.325	73	6.5
0.00	64	5.4

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 β Plate	Z0047311	2-8°C	30 Sep 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	M0430039	\leq -10°C	30 Sep 2022
Diluent 3	M0030097	\leq -10°C	31 Aug 2022
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
MSD Wash Buffer (20X), 100mL	Not Kit Specific
Plate Seals	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		27 MAY 2020