

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human MIP-1 $\beta$  Kit  
**Kit Catalog Numbers:** K151NR-Series; K156NR-Series  
**Kit Lot Number:** K0081165  
**Expiration Date:** 28 Feb 2021

### 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 $\beta$	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 $\beta$	1.93	80% - 120%	< 20%	750	80% - 120%	< 20%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	MIP-1 $\beta$	Calibrator	1240	pg/mL
		Chemokine Panel 1 (human) Control 1	991	pg/mL
		Chemokine Panel 1 (human) Control 2	172	pg/mL
		Chemokine Panel 1 (human) Control 3	25.5	pg/mL

Controls tested may not be the same lot as shipped.

### Standard Curve Data:

MIP-1 $\beta$ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1240	428,000	2.5
310	75,900	4.7
77.5	10,800	4.8
19.4	1,500	6.1
4.84	352	4.0
1.21	127	4.9
0.303	79	4.7
0.00	66	8.4

**CERTIFICATE OF ANALYSIS**

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1β Plate	Z0046992	2-8°C	30 Jun 2021
Anti-hu MIP-1β Antibody	D0081121	2-8°C	31 Jan 2022
Chemokine Panel 1 (human) Calibrator Blend	A0080183	2-8°C	31 Oct 2023
Diluent 43	M0430029	≤ -10°C	30 Apr 2021
Diluent 3	M0030085	≤ -10°C	28 Feb 2021
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

**Additional Comments:**

	Assay	Spot	Specification	Comment
1	MIP-1β	1	LLOQ CV	MIP-1β meets all specifications at LLOQ of 1.93 pg/mL.

**Statement:**

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		21 OCT 2019