

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human MIP-1 $\beta$   
**Kit Catalog Numbers:** K151NR-Series  
**Kit Lot Number:** K0081054  
**Expiration Date:** 31 Dec 2019

### 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	hu 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	hu MIP-1 $\beta$	1.02	80% - 120%	< 20%	750	80% - 120%	< 20%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu MIP-1 $\beta$	Calibrator	1180	pg/mL
		Chemokine Panel 1 (human) Control 1	1000	pg/mL
		Chemokine Panel 1 (human) Control 2	176	pg/mL
		Chemokine Panel 1 (human) Control 3	25.8	pg/mL

Controls tested may not be the same lot as shipped.

### Standard Curve Data:

hu MIP-1 $\beta$ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1180	476,000	4.0
295	93,700	2.7
73.8	16,000	4.1
18.4	2,510	2.9
4.61	534	4.1
1.15	178	4.4
0.288	105	4.8
0.00	75	8.7



## CERTIFICATE OF ANALYSIS

Page 2 of 2

### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 $\beta$ Plate	Z0046874	2-8°C	31 Dec 2020
Anti-hu MIP-1 $\beta$ Antibody	D0080991	2-8°C	31 Dec 2019
Chemokine Panel 1 (human) Calibrator Blend	A0080171	2-8°C	30 Jun 2023
Diluent 43	M0430027	$\leq$ -10°C	31 Dec 2020
Diluent 3	M0030082	$\leq$ -10°C	31 Oct 2020
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	Dawn Stollar	Quality Assurance		05 SEP 2018