

**CERTIFICATE OF ANALYSIS**

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human MIP-1β Kit  
**Kit Catalog Numbers:** K151NR-Series  
**Kit Lot Number:** K0080960  
**Expiration Date:** 31 Jan 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-1β	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-1β	2.27	80% - 120%	< 20%	750	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	hu MIP-1β	Calibrator	1190	pg/mL
		Chemokine Panel 1 (human) Control 1	940	pg/mL
		Chemokine Panel 1 (human) Control 2	158	pg/mL
		Chemokine Panel 1 (human) Control 3	26.9	pg/mL

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

hu MIP-1β - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1190	675,000	5.6
298	157,000	7.3
74.4	26,500	6.6
18.6	4,530	6.7
4.65	1,040	4.9
1.16	301	5.0
0.291	135	3.7
0.00	84	5.9

## CERTIFICATE OF ANALYSIS

Page 2 of 2

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 $\beta$ Plate	Z0046178	2-8°C	31 Jan 2019
Anti-hu MIP-1 $\beta$ Antibody	D0080861	2-8°C	30 Jun 2019
Chemokine Panel 1 (human) Calibrator Blend	A0080079	2-8°C	31 May 2021
Diluent 43	M0430021	$\leq$ -10°C	30 Sep 2019
Diluent 3	M0030077	$\leq$ -10°C	30 Nov 2019
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), 100 mL	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
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	02 APR 2018