

**CERTIFICATE OF ANALYSIS**

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human MIP-1 $\alpha$  Kit  
**Kit Catalog Numbers:** K151NQ-Series  
**Kit Lot Number:** K0081111  
**Expiration Date:** 31 Jan 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	hu MIP-1 $\alpha$	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 $\alpha$	13.8	80% - 120%	< 20%	743	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	hu MIP-1 $\alpha$	Calibrator	1060	pg/mL
		Chemokine Panel 1 (human) Control 1	929	pg/mL
		Chemokine Panel 1 (human) Control 2	411	pg/mL
		Chemokine Panel 1 (human) Control 3	167	pg/mL

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

hu MIP-1 $\alpha$ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1060	288,000	0.8
265	38,200	1.9
66.3	4,670	2.5
16.6	765	2.8
4.14	247	2.0
1.04	113	3.0
0.259	82	4.7
0.00	75	3.8

**CERTIFICATE OF ANALYSIS**

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 $\alpha$ Plate	Z0046941	2-8°C	31 Mar 2021
Anti-hu MIP-1 $\alpha$ Antibody	D0081123	2-8°C	31 Jan 2022
Chemokine Panel 1 (human) Calibrator Blend	A0080171	2-8°C	30 Jun 2023
Diluent 43	M0430028	$\leq$ -10°C	31 Jan 2021
Diluent 3	M0030085	$\leq$ -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)	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>	27 NOV 2018