



CERTIFICATE OF ANALYSIS

COA-6076

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human MIP-1 β Kit
Kit Catalog Numbers: K151NRD-1, K151NRD-2, K151NRD-4, K151NRG-1, K151NRG-2, K151NRG-4
Kit Lot Number: K0080384
Expiration Date: 31 May 2017

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 β	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 β	2.27	80% - 120%	< 20%	750	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu MIP-1 β	Calibrator	982	pg/mL
		Chemokine Panel 1 (human) Control 1	768	pg/mL
		Chemokine Panel 1 (human) Control 2	124	pg/mL
		Chemokine Panel 1 (human) Control 3	22.4	pg/mL

Controls tested may not be the same lot as shipped.

CERTIFICATE OF ANALYSIS

COA-6076

Page 2 of 2

Standard Curve Data:

hu MIP-1 β - Spot 1		
Conc. pg/mL	ECL Signal	%CV
982	361,000	4.6
246	73,800	4.3
61.4	11,800	5.6
15.3	2,530	4.9
3.84	606	4.3
0.959	205	3.9
0.240	96	4.1
0.00	63	7.2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 β Plate	Z0045808	2-8°C	31 Dec 2017
Anti-hu MIP-1 β Antibody	D0080534	2-8°C	31 Oct 2017
Chemokine Panel 1 (human) Calibrator Blend	A0080046	2-8°C	31 May 2017
Diluent 43	M0430014	$\leq -10^{\circ}\text{C}$	30 Sep 2017
Diluent 3	M0030067	$\leq -10^{\circ}\text{C}$	30 Nov 2017
MSD Read Buffer T (4X)	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 kit was tested using MSD procedures and it conforms to MSD's specifications. The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Matthew Chorley	Quality Assurance		30 OCT 2015