

CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-06935

Product Description: V-PLEX® and V-PLEX Plus Human MIP-1β Kit
Kit Catalog Numbers: K151NRD-1, K151NRD-2, K151NRD-4, K151NRG-1, K151NRG-2, K151NRG-4
Kit Lot Number: K0080563
Expiration Date: 31 Dec 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-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	1010	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

Page 2 of 2

COA-06935

Standard Curve Data:

hu MIP-1 β - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1010	439,000	2.1
253	89,500	5.6
63.1	15,200	2.9
15.8	2,960	2.6
3.95	691	1.8
0.986	225	2.5
0.247	108	3.4
0.00	72	6.3


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 β Plate	Z0046116	2-8°C	30 Sep 2018
Chemokine Panel 1 (human) Calibrator Blend	A0080067	2-8°C	31 Jan 2018
Diluent 43	M0430015	\leq -10°C	31 Dec 2017
Diluent 3	M0030069	\leq -10°C	30 Apr 2018
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), 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 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	Dawn Stollar	Quality Assurance		31 MAY 2016