

CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-07326

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: K0080592
Expiration Date: 31 Jan 2018

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-07326

Standard Curve Data:

hu MIP-1 β - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1010	528,000	2.7
253	110,000	3.5
63.1	19,900	3.4
15.8	3,990	2.3
3.95	969	4.0
0.986	291	3.7
0.247	129	4.6
0.00	74	5.5


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 β Plate	Z0046178	2-8°C	31 Jan 2019
Anti-hu MIP-1 β Antibody	D0080685	2-8°C	30 Nov 2018
Chemokine Panel 1 (human) Calibrator Blend	A0080067	2-8°C	31 Jan 2018
Diluent 43	M0430017	\leq -10°C	31 Dec 2018
Diluent 3	M0030070	\leq -10°C	31 Jul 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		24 SEP 2016