

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human MIP-1 β Kit
Kit Catalog Numbers: K151NR-Series
Kit Lot Number: K0080989
Expiration Date: 30 Sep 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-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	1.02	80% - 120%	< 20%	750	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu MIP-1 β	Calibrator	1110	pg/mL
		Chemokine Panel 1 (human) Control 1	1020	pg/mL
		Chemokine Panel 1 (human) Control 2	175	pg/mL
		Chemokine Panel 1 (human) Control 3	25.9	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu MIP-1 β - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1110	418,000	6.2
278	90,800	4.1
69.4	15,900	2.2
17.3	2,710	2.2
4.34	627	3.6
1.08	195	3.2
0.271	100	2.9
0.00	69	4.1

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 β Plate	Z0046799	2-8°C	30 Sep 2020
Anti-hu MIP-1 β Antibody	D0080898	2-8°C	30 Sep 2019
Chemokine Panel 1(Human) Calibrator Blend	A0080122	2-8°C	28 Feb 2022
Diluent 43	M0430021	\leq -10°C	30 Sep 2019
Diluent 3	M0030076	\leq -10°C	30 Sep 2019
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Wash Buffer (20X), 100mL	Not Kit Specific
Plate Seals	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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	04 JUN 2018