

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-Plex Plus Human MIP-1β Kit
Kit Catalog Numbers: K151NR-Series
Kit Lot Number: K0081123
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-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	1.02	80% - 120%	< 20%	750	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu MIP-1β	Calibrator	1180	pg/mL
		Chemokine Panel 1 (human) Control 1	1000	pg/mL
		Chemokine Panel 1 (human) Control 2	176	pg/mL
		Chemokine Panel 1 (human) Control 3	25.8	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu MIP-1β - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1180	386,000	6.2
295	64,400	7.8
73.8	8,820	8.6
18.4	1,190	8.8
4.61	247	5.9
1.15	106	5.4
0.288	80	6.6
0.00	71	4.2

CERTIFICATE OF ANALYSIS

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 β Plate	Z0046944	2-8°C	31 Mar 2021
Anti-hu MIP-1 β Antibody	D0081121	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), 100 mL	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>	28 NOV 2018