

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human MDC Kit
Kit Catalog Numbers: K151NP-Series
Kit Lot Number: K0080996
Expiration Date: 30 Jun 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 MDC	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu MDC	88.3	80% - 120%	< 20%	7500	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu MDC	Calibrator	11600	pg/mL
		Chemokine Panel 1 (human) Control 1	9420	pg/mL
		Chemokine Panel 1 (human) Control 2	2770	pg/mL
		Chemokine Panel 1 (human) Control 3	878	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu MDC - Spot 1		
Conc. pg/mL	ECL Signal	%CV
11600	740,000	3.8
2900	103,000	3.3
725	16,200	3.5
181	3,320	2.5
45.3	863	2.1
11.3	257	3.5
2.83	116	3.2
0.00	64	3.8

CERTIFICATE OF ANALYSIS

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MDC Plate	Z0046803	2-8°C	30 Sep 2020
Anti-hu MDC Antibody	D0080866	2-8°C	30 Jun 2019
Chemokine Panel 1 (human) Calibrator Blend	A0080159	2-8°C	31 Dec 2022
Diluent 43	M0430024	≤ -10°C	31 Jan 2020
Diluent 3	M0030081	≤ -10°C	30 Sep 2020
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>	23 MAY 2018