

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human MIP-1 α Kit
Kit Catalog Numbers: K151NQ-Series
Kit Lot Number: K0080967
Expiration Date: 31 Jul 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 α	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 α	13.8	80% - 120%	< 20%	743	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu MIP-1 α	Calibrator	1020	pg/mL
		Chemokine Panel 1 (human) Control 1	896	pg/mL
		Chemokine Panel 1 (human) Control 2	409	pg/mL
		Chemokine Panel 1 (human) Control 3	151	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu MIP-1 α - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1020	231,000	2.6
255	27,600	2.3
63.8	2,390	3.7
15.9	262	4.6
3.98	85	4.6
0.996	68	5.1
0.249	62	5.2
0.00	62	4.2

CERTIFICATE OF ANALYSIS

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 α Plate	Z0046454	2-8°C	30 Sep 2019
Anti-hu MIP-1 α Antibody	D0080902	2-8°C	31 Jul 2019
Chemokine Panel 1 (human) Calibrator Blend	A0080159	2-8°C	31 Dec 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

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>	13 JUN 2018