

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, K156NQ-Series
Kit Lot Number: K0081203
Expiration Date: 30 Jun 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	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	MIP-1 α	13.8	80% - 120%	< 20%	743	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	MIP-1 α	Calibrator	1090	pg/mL
		Chemokine Panel 1 (human) Control 1	920	pg/mL
		Chemokine Panel 1 (human) Control 2	403	pg/mL
		Chemokine Panel 1 (human) Control 3	169	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

MIP-1 α - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1090	216,000	1.6
273	25,100	3.2
68.1	2,110	4.0
17.0	228	3.2
4.26	82	4.5
1.06	68	9.0
0.266	64	3.4
0.00	67	5.5

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 α Plate	Z0047040	2-8°C	31 Aug 2021
Anti-hu MIP-1 α Antibody	D0081123	2-8°C	31 Jan 2022
Chemokine Panel 1 (human) Calibrator Blend	A0080183	2-8°C	31 Oct 2023
Diluent 43	M0430030	≤ -10°C	30 Jun 2021
Diluent 3	M0030087	≤ -10°C	30 Jun 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), 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 product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality		17 APR 2019