

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Mouse IP-10 Kit  
**Kit Catalog Numbers:** K152NV-Series  
**Kit Lot Number:** K0081156  
**Expiration Date:** 31 Jan 2021

### Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Cytokine Panel 1 (mouse) Control 1		Cytokine Panel 1 (mouse) Control 2		Cytokine Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	ms IP-10	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	ms IP-10	2.15	75% - 125%	< 20%	650	75% - 125%	< 20%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms IP-10	Calibrator	910	pg/mL
		Cytokine Panel 1 (mouse) Control 1	947	pg/mL
		Cytokine Panel 1 (mouse) Control 2	245	pg/mL
		Cytokine Panel 1 (mouse) Control 3	77.9	pg/mL

Controls tested may not be the same lot as shipped.

### Standard Curve Data:

ms IP-10 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
910	340,000	4.4
228	83,500	4.7
56.9	20,900	3.8
14.2	5,210	4.7
3.55	1,380	2.4
0.889	419	2.2
0.222	185	2.7
0.00	111	4.7

## CERTIFICATE OF ANALYSIS

Page 2 of 2


### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IP-10 Plate	Z0046978	2-8°C	30 Apr 2021
Anti-ms IP-10 Antibody	D0081117	2-8°C	31 Jan 2022
Cytokine Panel 1 (mouse) Calibrator Blend	A0080135	2-8°C	30 Jun 2022
Diluent 41	M0410025	≤ -10°C	31 Jan 2021
Diluent 45	M0450023	≤ -10°C	31 Jan 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
Cytokine Panel 1 (mouse) Control 1	Not Kit Specific
Cytokine Panel 1 (mouse) Control 2	Not Kit Specific
Cytokine Panel 1 (mouse) 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
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		21 DEC 2018