

**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:** K0081068

**Expiration Date:** 30 Nov 2019

**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	235,000	2.5
228	56,700	2.8
56.9	13,400	3.5
14.2	3,360	3.4
3.55	944	1.4
0.889	420	24.0
0.222	212	4.8
0.00	155	1.8

## CERTIFICATE OF ANALYSIS

Page 2 of 2

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IP-10 Plate	Z0046889	2-8°C	31 Jan 2021
Anti-ms IP-10 Antibody	D0080977	2-8°C	30 Nov 2019
Cytokine Panel 1 (mouse) Calibrator Blend	A0080135	2-8°C	30 Jun 2022
Diluent 41	M0410024	≤ -10°C	30 Sep 2020
Diluent 45	M0450021	≤ -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
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	<i>Dawn Stollar</i>	12 SEP 2018