

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: K0081083
Expiration Date: 30 Sep 2020

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	75% - 125%	< 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	288,000	3.0
228	72,300	2.9
56.9	17,400	3.3
14.2	4,600	2.8
3.55	1,220	3.4
0.889	418	3.4
0.222	211	4.9
0.00	143	2.8

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IP-10 Plate	Z0046907	2-8°C	28 Feb 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	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), 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
Review/Approval	Dawn Stollar	Quality Assurance		08 OCT 2018