

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: K0080936
Expiration Date: 31 Aug 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	883	pg/mL
		Cytokine Panel 1 (mouse) Control 2	251	pg/mL
		Cytokine Panel 1 (mouse) Control 3	82.7	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	351,000	2.5
228	94,600	1.7
56.9	23,600	2.1
14.2	5,930	2.1
3.55	1,590	3.5
0.889	482	3.5
0.222	208	2.9
0.00	116	2.4

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IP-10 Plate	Z0046716	2-8°C	30 Jun 2020
Anti-ms IP-10 Antibody	D0080939	2-8°C	31 Aug 2019
Cytokine Panel 1 (mouse) Calibrator Blend	A0080135	2-8°C	30 Jun 2022
Diluent 41	M0410020	≤ -10°C	28 Feb 2020
Diluent 45	M0450019	≤ -10°C	31 Dec 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
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		27 FEB 2018