

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: K0081343
Expiration Date: 31 Dec 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	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	IP-10	2.15	75% - 125%	< 20%	650	75% - 125%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IP-10	Calibrator	1100	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:

IP-10 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1100	265,000	4.6
275	65,300	3.4
68.8	15,600	3.5
17.2	3,960	4.7
4.30	1,120	4.2
1.07	369	3.6
0.269	192	3.9
0.00	132	3.9

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IP-10 Plate	Z0047197	2-8°C	31 Mar 2022
Anti-ms IP-10 Antibody	D0081383	2-8°C	31 May 2023
Cytokine Panel 1 (mouse) Calibrator Blend	A0080218	2-8°C	31 Oct 2024
Diluent 41	M0410029	≤ -10°C	31 Dec 2021
Diluent 45	M0450030	≤ -10°C	28 Feb 2022
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	Karen Hamilla	Quality		21 JAN 2020