

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Mouse IP-10 Kit  
**Kit Lot Number:** K0082292  
**Expiration Date:** 30 Jun 2025  
**Kit Catalog Numbers:** K152NV-Series

**Precision and Accuracy 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 Calibrator:**


Spot	Name	Conc.	Units
1	IP-10	1010	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IP-10 SECTOR Plate	Z0047864	2-8°C	31 Oct 2025
Anti-ms IP-10 Antibody	D0081937	2-8°C	31 Mar 2026
Cytokine Panel 1 (mouse) Calibrator Blend	A0080353	2-8°C	29 Feb 2028
Diluent 41	M0410052	≤ -10°C	30 Sep 2025
Diluent 45	M0450051	≤ -10°C	30 Jun 2025

Additional components may be provided to execute the assay. Additional components are not lot-specific.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		26 JUL 2023