

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX® and V-PLEX Plus Mouse IP-10 Kit**Lot Number:** K0082634**Expiration Date:** 30 NOV 2026**Catalog Number:** K152NV-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-ms IP-10 Antibody	D0082136	2-8°C	31 JAN 2027
Cytokine Panel 1 (mouse) Calibrator Blend	A0080353	2-8°C	29 FEB 2028
Mouse IP-10, 96-Well Small Spot SECTOR Plate	Z0048068	2-8°C	31 JUL 2027
Diluent 41	M0410056	≤ -10°C	31 MAR 2027
Diluent 45	M0450054	≤ -10°C	30 NOV 2026

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

**Precision, Accuracy, and Sensitivity 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

Dilution to top of Curve: 1X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		05 MAR 2025