



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Mouse IFN- γ Kit**Lot Number:** K0082821**Expiration Date:** 30 NOV 2027**Catalog Number:** K152QO-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-ms IFN- γ Antibody	D0082263	2-8°C	30 NOV 2027
Mouse IFN- γ 96-Well Small Spot SECTOR Plate	Z0048193	2-8°C	30 APR 2028
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080440	2-8°C	31 AUG 2030
Diluent 41	M0410059	$\leq -10^{\circ}\text{C}$	31 DEC 2027
Diluent 45	M0450056	$\leq -10^{\circ}\text{C}$	29 FEB 2028

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

Precision, Accuracy, and Sensitivity Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (mouse) Control 1		Proinflammatory Panel 1 (mouse) Control 2		Proinflammatory Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IFN- γ	80% - 120%	< 20%	90% - 140%	< 25%	90% - 140%	< 25%	90% - 140%	< 25%


Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IFN- γ	0.39	75% - 125%	< 25%	570	75% - 125%	< 25%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IFN- γ	991	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		23 DEC 2025