



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Mouse IFN- γ Kit**Lot Number:** K0082963**Expiration Date:** 31 DEC 2027**Catalog Number:** K152QO-Series**Lot-Specific Components:**

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Anti-ms IFN- γ Antibody	D0082443	2-8°C	31 AUG 2028
Mouse IFN- γ 96-Well Small Spot SECTOR Plate	Z0048270	2-8°C	30 NOV 2028
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080429	2-8°C	31 MAY 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:

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

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	IFN- γ	0.390	75% - 125%	< 25%	570	75% - 125%	< 25%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IFN- γ	938	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		28 JUL 2026