



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Human IFN- γ Kit**Lot Number:** K0082692**Expiration Date:** 30 SEP 2026**Catalog Number:** K151QO-Series, K156QO-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IFN- γ Antibody	D0082411	2-8°C	31 JUL 2028
Human IFN- γ Plate	Z0048111	2-8°C	30 SEP 2027
Proinflammatory Panel 1 (human) Calibrator Blend	A0080427	2-8°C	28 FEB 2030
Diluent 2	M0020130	$\leq -10^{\circ}\text{C}$	30 SEP 2026
Diluent 3	M0030140	$\leq -10^{\circ}\text{C}$	30 SEP 2027

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

Precision, Accuracy, and Sensitivity Specifications:

		Calibrator		Proinflammatory Panel 1 (human) Control 1		Proinflammatory Panel 1 (human) Control 2		Proinflammatory Panel 1 (human) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IFN- γ	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	IFN- γ	1.76	80% - 120%	< 20%	938	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IFN- γ	1420	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		08 MAY 2025