



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Human IFN- γ Kit, QuickPlex
Lot Number: K0082646
Expiration Date: 31 DEC 2026
Catalog Number: K151QO-Series; K156QO-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IFN- γ Antibody	D0082162	2-8°C	30 APR 2027
Human IFN- γ 96-Well Small Spot QuickPlex Ultra Plate	Z0048076	2-8°C	31 JUL 2027
Proinflammatory Panel 1 (human) Calibrator Blend	A0080412	2-8°C	31 DEC 2029
Diluent 2	M0020131	$\leq -10^{\circ}\text{C}$	31 DEC 2026
Diluent 3	M0030136	$\leq -10^{\circ}\text{C}$	31 MAY 2027

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

Spot	Name	LLOQ			ULOQ		
		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	1440	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		18 MAR 2025