



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Human IFN- γ Kit
Lot Number: K0082782
Expiration Date: 30 NOV 2027
Catalog Number: K151QO-Series, K156QO-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IFN- γ Antibody	D0082440	2-8°C	31 AUG 2028
Human IFN- γ 96-Well Small Spot SECTOR Plate	Z0048169	2-8°C	29 FEB 2028
Proinflammatory Panel 1 (human) Calibrator Blend	A0080436	2-8°C	30 JUN 2030
Diluent 2	M0020135	$\leq -10^{\circ}\text{C}$	30 NOV 2027
Diluent 3	M0030141	$\leq -10^{\circ}\text{C}$	30 NOV 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	1290	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		02 OCT 2025