



### CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX® and V-PLEX Plus Human IFN- $\gamma$  Kit  
**Lot Number:** K0082542  
**Expiration Date:** 31 MAR 2026  
**Catalog Number:** K151QO-Series, K156QO-Series

#### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IFN- $\gamma$ Antibody	D0082286	2-8°C	31 DEC 2027
Human IFN- $\gamma$ Plate	Z0048019	2-8°C	31 JAN 2027
Proinflammatory Panel 1 (human) Calibrator Blend	A0080400	2-8°C	30 JUN 2029
Diluent 2	M0020128	$\leq -10^\circ\text{C}$	31 MAR 2026
Diluent 3	M0030129	$\leq -10^\circ\text{C}$	30 NOV 2026

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- $\gamma$	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%

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

#### Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IFN- $\gamma$	1060	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
<b>Review/Approval</b>	Karen Hamilla	Quality		01 OCT 2024



**CERTIFICATE OF ANALYSIS**