

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX® and V-PLEX Plus Human IFN- $\gamma$  Kit**Lot Number:** K0082622**Expiration Date:** 31 MAR 2027**Catalog Number:** K151QO-Series, K156QO-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IFN- $\gamma$ Antibody	D0082356	2-8°C	31 MAR 2028
Human IFN- $\gamma$ 96-Well Small Spot SECTOR Plate	Z0048060	2-8°C	30 JUN 2027
Proinflammatory Panel 1 (human) Calibrator Blend	A0080385	2-8°C	31 DEC 2028
Diluent 2	M0020132	$\leq -10^\circ\text{C}$	31 MAR 2027
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- $\gamma$	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- $\gamma$	1380	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		07 APR 2025