

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human IFN- $\gamma$  Kit  
**Kit Lot Number:** K0082214  
**Expiration Date:** 31 Jul 2024  
**Kit Catalog Numbers:** K151QO-Series, K156QO-Series

**Precision and Accuracy 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%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 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- $\gamma$	1470	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IFN- $\gamma$ Plate	Z0047807	2-8°C	31 Jul 2025
Anti-hu IFN- $\gamma$ Antibody	D0081918	2-8°C	30 Apr 2026
Proinflammatory Panel 1 (human) Calibrator Blend	A0080345	2-8°C	31 Oct 2027
Diluent 2	M0020124	$\leq$ -10°C	28 Feb 2025
Diluent 3	M0030113	$\leq$ -10°C	31 Jul 2024

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

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		08 MAR 2023