

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-Plex® and V-Plex Plus Human IFN- $\gamma$  Kit  
**Kit Lot Number:** K0081701  
**Expiration Date:** 31 May 2023  
**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- $\gamma$	1530	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IFN- $\gamma$ Plate	Z0047521	2-8°C	30 Sep 2023
Anti-hu IFN- $\gamma$ Antibody	D0081642	2-8°C	30 Sep 2024
Proinflammatory Panel 1 (human) Calibrator Blend	A0080289	2-8°C	31 Mar 2026
Diluent 2	M0020115	$\leq$ -10°C	31 May 2023
Diluent 3	M0030107	$\leq$ -10°C	30 Sep 2023

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		14 JUN 2021