

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IFN- γ Kit
Kit Lot Number: K0082153
Expiration Date: 28 Feb 2025
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- γ	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 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	1520	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IFN- γ Plate	Z0047778	2-8°C	31 Mar 2025
Anti-hu IFN- γ Antibody	D0081875	2-8°C	28 Feb 2026
Proinflammatory Panel 1 (human) Calibrator Blend	A0080334	2-8°C	31 Jul 2027
Diluent 2	M0020123	\leq -10°C	28 Feb 2025
Diluent 3	M0030118	\leq -10°C	31 Mar 2025

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 NOV 2022