

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IFN- γ Kit

Kit Lot Number: K0082415

Expiration Date: 30 Jun 2026

Kit Catalog Numbers: K151QO-Series; K156QO-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IFN- γ Plate	Z0047947	2-8°C	30 Jun 2026
Anti-hu IFN- γ Antibody	D0082162	2-8°C	30 Apr 2027
Proinflammatory Panel 1 (human) Calibrator Blend	A0080382	2-8°C	30 Nov 2028
Diluent 2	M0020129	$\leq -10^\circ\text{C}$	30 Jun 2026
Diluent 3	M0030126	$\leq -10^\circ\text{C}$	31 Jul 2026

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

Precision and Accuracy Specifications:

		Calibrator		Proinflammatory Panel 1 (human) Control 1		Proinflammatory Panel 1 (human) Control 2		Proinflammatory Panel 1 (human) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IFN- γ	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%


		LLOQ			ULOQ		
Spot	Name	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	1340	pg/mL

Dilution to Top of Curve = 1X

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

	Name	Function	Signature	Date
Review/Approval	Michelle Likiam	Quality		26 Mar 2024