

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IFN- γ Kit
Kit Lot Number: K0081578
Expiration Date: 31 Dec 2022
Kit Catalog Numbers: K151QO-Series, K156QO-Series

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%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 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	1360	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IFN- γ Plate	Z0047409	2-8°C	31 Mar 2023
Anti-hu IFN- γ Antibody	D0081528	2-8°C	29 Feb 2024
Proinflammatory Panel 1 (human) Calibrator Blend	A0080267	2-8°C	31 Aug 2025
Diluent 2	M0020114	$\leq -10^{\circ}\text{C}$	31 Mar 2023
Diluent 3	M0030101	$\leq -10^{\circ}\text{C}$	31 Dec 2022

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	Xiomara Talavera	Quality		01 DEC 2020