

CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Kit Lot Number: K0082090

Expiration Date: 31 Dec 2023

Kit Catalog Numbers: K152QO-Series

Precision and Accuracy Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (mouse) Control 1		Proinflammatory Panel 1 (mouse) Control 2		Proinflammatory Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IFN- γ	80% - 120%	< 20%	90% - 140%	< 25%	90% - 140%	< 25%	90% - 140%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IFN- γ	0.390	75% - 125%	< 25%	570	75% - 125%	< 25%

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	IFN- γ	742	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IFN- γ Plate	Z0047742	2-8°C	31 Jan 2025
Anti-ms IFN- γ Antibody	D0081660	2-8°C	31 Oct 2024
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080282	2-8°C	31 Mar 2026
Diluent 41	M0410044	\leq -10°C	30 Apr 2024
Diluent 45	M0450042	\leq -10°C	31 Dec 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		25 AUG 2022