

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Mouse IFN- $\gamma$  Kit  
**Kit Lot Number:** K0082391  
**Expiration Date:** 30 Nov 2025  
**Kit Catalog Numbers:** K152QO-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IFN- $\gamma$ Plate	Z0047926	2-8°C	30 Apr 2026
Anti-ms IFN- $\gamma$ Antibody	D0081921	2-8°C	28 Feb 2026
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080351	2-8°C	31 Dec 2027
Diluent 41	M0410053	$\leq$ -10°C	30 Nov 2025
Diluent 45	M0450053	$\leq$ -10°C	31 Mar 2026

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

**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- $\gamma$	80% - 120%	< 20%	90% - 140%	< 25%	90% - 140%	< 25%	90% - 140%	< 25%


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

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	IFN- $\gamma$	1030	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		07 Dec 2023