

**CERTIFICATE OF ANALYSIS**

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Mouse IFN- $\gamma$  Kit

**Kit Catalog Numbers:** K152QO-Series

**Kit Lot Number:** K0081542

**Expiration Date:** 31 Dec 2022

**Acceptance Data/ Percent Recovery 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 Material Tested :**

Spot	Name	Item	Conc	Units
1	IFN- $\gamma$	Calibrator	785	pg/ml
		Proinflammatory Panel 1 (mouse) Control 1	469	pg/ml
		Proinflammatory Panel 1 (mouse) Control 2	32	pg/ml
		Proinflammatory Panel 1 (mouse) Control 3	2.12	pg/ml

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

IFN- $\gamma$ - Spot 1		
Conc. pg/ml	ECL Signal	%CV
785	348,000	2.0
196	82,500	1.7
49.1	20,200	1.1
12.3	5,010	2.0
3.07	1,280	2.5
0.767	360	1.9
0.192	138	3.2
0.00	64	6.4

## CERTIFICATE OF ANALYSIS

Page 2 of 2


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IFN- $\gamma$ Plate	Z0047384	2-8°C	28 Feb 2023
Anti-ms IFN- $\gamma$ Antibody	D0081487	2-8°C	31 Oct 2023
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080252	2-8°C	30 Apr 2025
Diluent 41	M0410034	$\leq -10^{\circ}\text{C}$	31 Dec 2022
Diluent 45	M0450035	$\leq -10^{\circ}\text{C}$	31 Jan 2023
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Plate Seals	Not Kit Specific
Wash Buffer (20X), 100mL	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 1	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 2	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 3	Not Kit Specific

**Statement:**

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		08 OCT 2020