

**CERTIFICATE OF ANALYSIS**

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human IFN- $\gamma$  Kit  
**Kit Catalog Numbers:** K151QO-Series, K156QO-Series  
**Kit Lot Number:** K0081443  
**Expiration Date:** 28 Feb 2022

**Acceptance Data/ Percent Recovery Specifications:**

Spot	Name	Calibrator		Proinflammatory Panel 1 (human) Control 1		Proinflammatory Panel 1 (human) Control 2		Proinflammatory Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IFN- $\gamma$	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IFN- $\gamma$	1.76	80% - 120%	< 20%	938	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	IFN- $\gamma$	Calibrator	1470	pg/mL
		Proinflammatory Panel 1 (human) Control 1	530	pg/mL
		Proinflammatory Panel 1 (human) Control 2	145	pg/mL
		Proinflammatory Panel 1 (human) Control 3	35.3	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
1470	74,400	1.5
368	18,600	1.7
91.9	4,580	2.3
23.0	1,180	1.5
5.74	339	3.2
1.44	131	2.8
0.359	78	6.8
0.00	65	7.9

**CERTIFICATE OF ANALYSIS**


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IFN- $\gamma$ Plate	Z0047283	2-8°C	31 Aug 2022
Anti-hu IFN- $\gamma$ Antibody	D0081180	2-8°C	31 Mar 2022
Proinflammatory Panel 1 (human) Calibrator Blend	A0080229	2-8°C	30 Sep 2024
Diluent 2	M0020109	$\leq -10^{\circ}\text{C}$	28 Feb 2022
Diluent 3	M0030096	$\leq -10^{\circ}\text{C}$	30 Jun 2022
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), 100 mL	Not Kit Specific
Proinflammatory Panel 1 (human) Control 1	Not Kit Specific
Proinflammatory Panel 1 (human) Control 2	Not Kit Specific
Proinflammatory Panel 1 (human) 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		01 JUL 2020