

## 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:** K0081298  
**Expiration Date:** 31 Dec 2021

### 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	1280	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
1280	107,000	2.3
320	26,900	2.9
80.0	6,670	3.4
20.0	1,700	3.2
5.00	474	3.5
1.25	175	4.2
0.313	95	2.6
0.00	71	6.8

**CERTIFICATE OF ANALYSIS**


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IFN- $\gamma$ Plate	Z0047151	2-8°C	31 Jan 2022
Anti-hu IFN- $\gamma$ Antibody	D0081304	2-8°C	31 Oct 2022
Proinflammatory Panel 1 (human) Calibrator Blend	A0080216	2-8°C	30 Jun 2024
Diluent 2	M0020109	$\leq$ -10°C	28 Feb 2022
Diluent 3	M0030090	$\leq$ -10°C	31 Dec 2021
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 (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>	Xiomara Talavera	Quality		09 DEC 2019