

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human IFN- $\gamma$  Kit  
**Kit Catalog Numbers:** K151QO-Series  
**Kit Lot Number:** K0081093  
**Expiration Date:** 31 Jan 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	hu 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	hu IFN- $\gamma$	1.76	80% - 120%	< 20%	938	80% - 120%	< 20%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IFN- $\gamma$	Calibrator	1250	pg/mL
		Proinflammatory Panel 1 (human) Control 1	512	pg/mL
		Proinflammatory Panel 1 (human) Control 2	132	pg/mL
		Proinflammatory Panel 1 (human) Control 3	40.9	pg/mL

Controls tested may not be the same lot as shipped.

### Standard Curve Data:

hu IFN- $\gamma$ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1250	71,300	2.1
313	17,100	6.6
78.1	4,290	3.4
19.5	1,110	2.9
4.88	326	3.8
1.22	134	3.1
0.305	89	6.1
0.00	73	3.5

## CERTIFICATE OF ANALYSIS


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IFN- $\gamma$ Plate	Z0046918	2-8°C	28 Feb 2021
Anti-hu IFN- $\gamma$ Antibody	D0081074	2-8°C	31 Aug 2021
Proinflammatory Panel 1 (human) Calibrator Blend	A0080180	2-8°C	30 Sep 2023
Diluent 2	M0020102	$\leq$ -10°C	28 Feb 2021
Diluent 3	M0030084	$\leq$ -10°C	31 Jan 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), 1000mL	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>	Dawn Stollar	Quality Assurance		16 JAN 2019