

**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:** K0080977

**Expiration Date:** 30 Sep 2019

**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- $\gamma$	7.47	80% - 120%	< 20%	938	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	hu IFN- $\gamma$	Calibrator	1460	pg/mL
		Proinflammatory Panel 1 (human) Control 1	682	pg/mL
		Proinflammatory Panel 1 (human) Control 2	155	pg/mL
		Proinflammatory Panel 1 (human) Control 3	52.3	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
1460	102,000	1.4
365	24,800	2.6
91.3	6,350	2.4
22.8	1,620	1.3
5.70	464	2.3
1.43	174	2.4
0.356	104	3.0
0.00	79	4.3

## CERTIFICATE OF ANALYSIS

Page 2 of 2

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IFN- $\gamma$ Plate	Z0046785	2-8°C	31 Aug 2020
Anti-hu IFN- $\gamma$ Antibody	D0081074	2-8°C	31 Aug 2021
Proinflammatory Panel 1 (human) Calibrator Blend	A0080153	2-8°C	31 Oct 2022
Diluent 2	M0020098	$\leq$ -10°C	28 Feb 2020
Diluent 3	M0030076	$\leq$ -10°C	30 Sep 2019
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>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	24 APR 2018