

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IFN-γ Kit

Kit Catalog Numbers: K151QO-Series

Kit Lot Number: K0080976

Expiration Date: 31 Jan 2020

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-γ	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-γ	7.47	80% - 120%	< 20%	938	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IFN-γ	Calibrator	1460	pg/mL
		Proinflammatory Panel 1 (human) Control 1	682	pg/mL
		Proinflammatory Panel 1 (human) Control 2	164	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-γ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1460	98,000	1.7
365	24,900	2.2
91.3	6,150	5.5
22.8	1,570	5.0
5.70	452	1.7
1.43	162	3.2
0.356	90	4.5
0.00	69	12.0

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IFN- γ Plate	Z0046650	2-8°C	28 Feb 2020
Anti-hu IFN- γ Antibody	D0081014	2-8°C	31 Jan 2020
Proinflammatory Panel 1 (human) Calibrator Blend	A0080139	2-8°C	31 Jul 2022
Diluent 2	M0020098	$\leq -10^{\circ}\text{C}$	28 Feb 2020
Diluent 3	M0030078	$\leq -10^{\circ}\text{C}$	31 Jan 2020
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)	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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	11 SEP 2018