

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IFN- γ Kit
Kit Catalog Numbers: K151QO-Series, K156QO-Series
Kit Lot Number: K0081400
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- γ	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

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

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IFN- γ	Calibrator	1470	pg/mL
		Proinflammatory Panel 1 (human) Control 1	568	pg/mL
		Proinflammatory Panel 1 (human) Control 2	173	pg/mL
		Proinflammatory Panel 1 (human) Control 3	36	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IFN- γ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1470	132,000	4.2
368	33,300	3.2
91.9	8,250	3.4
23.0	2,120	4.4
5.74	579	3.2
1.44	199	2.2
0.359	110	4.6
0.00	85	5.9

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IFN- γ Plate	Z0047242	2-8°C	30 Jun 2022
Anti-hu IFN- γ Antibody	D0081379	2-8°C	31 Mar 2023
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	M0030094	$\leq -10^{\circ}\text{C}$	31 May 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), 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
Review/Approval	Karen Hamilla	Quality		12 FEB 2020