

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: K0081389

Expiration Date: 30 Sep 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- 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	1190	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- γ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1190	100,000	1.6
298	25,300	1.5
74.4	6,290	1.8
18.6	1,610	1.5
4.65	466	2.1
1.16	168	0.8
0.291	99	6.7
0.00	75	8.2

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IFN- γ Plate	Z0047070	2-8°C	30 Sep 2021
Anti-hu IFN- γ Antibody	D0081254	2-8°C	31 Jul 2022
Proinflammatory Panel 1 (human) Calibrator Blend	A0080205	2-8°C	31 Mar 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
Review/Approval	Karen Hamilla	Quality		20 DEC 2019