

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: K0081516

Expiration Date: 31 Oct 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	1740	pg/mL
		Proinflammatory Panel 1 (human) Control 1	612	pg/mL
		Proinflammatory Panel 1 (human) Control 2	164	pg/mL
		Proinflammatory Panel 1 (human) Control 3	44.2	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
1740	131,000	2.7
435	33,200	3.2
109	8,470	2.6
27.2	2,170	2.5
6.80	590	2.2
1.70	198	2.1
0.425	102	5.2
0.00	72	6.6

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IFN- γ Plate	Z0047364	2-8°C	31 Dec 2022
Anti-hu IFN- γ Antibody	D0081474	2-8°C	30 Nov 2023
Proinflammatory Panel 1 (human) Calibrator Blend	A0080265	2-8°C	31 Jul 2025
Diluent 2	M0020113	\leq -10°C	30 Nov 2022
Diluent 3	M0030100	\leq -10°C	31 Oct 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		15 SEP 2020