

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-2 Kit  
**Kit Catalog Numbers:** K151QQ - Series, K156QQ - Series  
**Kit Lot Number:** K0081384  
**Expiration Date:** 31 Jan 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	IL-2	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-2	0.890	80% - 120%	< 20%	938	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	IL-2	Calibrator	1500	pg/mL
		Proinflammatory Panel 1 (human) Control 1	419	pg/mL
		Proinflammatory Panel 1 (human) Control 2	39.6	pg/mL
		Proinflammatory Panel 1 (human) Control 3	3.66	pg/mL

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

IL-2 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1500	453,000	1.8
375	129,000	2.0
93.8	34,300	2.1
23.4	8,890	2.2
5.86	2,310	4.8
1.46	623	2.7
0.366	220	3.6
0.00	84	3.5

## CERTIFICATE OF ANALYSIS

Page 2 of 2


### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-2 Plate	Z0047236	2-8°C	31 May 2022
Anti-hu IL-2 Antibody	D0081129	2-8°C	31 Jan 2022
Proinflammatory Panel 1 (human) Calibrator Blend	A0080216	2-8°C	30 Jun 2024
Diluent 2	M0020109	≤ -10°C	28 Feb 2022
Diluent 3	M0030094	≤ -10°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
<b>Review/Approval</b>	Karen Hamilla	Quality		18 FEB 2020