

## 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:** K0081394  
**Expiration Date:** 30 Sep 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	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	1420	pg/mL
		Proinflammatory Panel 1 (human) Control 1	409	pg/mL
		Proinflammatory Panel 1 (human) Control 2	39.4	pg/mL
		Proinflammatory Panel 1 (human) Control 3	3.49	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
1420	432,000	2.3
355	119,000	3.0
88.8	31,500	2.5
22.2	7,960	2.3
5.55	2,110	2.0
1.39	586	3.7
0.347	215	3.0
0.00	91	4.8

**CERTIFICATE OF ANALYSIS**


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-2 Plate	Z0046807	2-8°C	30 Sep 2020
Anti-hu IL-2 Antibody	D0081096	2-8°C	30 Nov 2021
Proinflammatory Panel 1 (human) Calibrator Blend	A0080158	2-8°C	31 May 2023
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		10 MAR 2020