

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-2 Kit  
**Kit Catalog Numbers:** K151QQ-Series  
**Kit Lot Number:** K0080905  
**Expiration Date:** 31 Jan 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	hu 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	hu IL-2	0.890	80% - 120%	< 20%	938	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	hu IL-2	Calibrator	1520	pg/mL
		Proinflammatory Panel 1 (human) Control 1	431	pg/mL
		Proinflammatory Panel 1 (human) Control 2	42.3	pg/mL
		Proinflammatory Panel 1 (human) Control 3	3.83	pg/mL

Controls tested may not be the same lot as shipped.

## CERTIFICATE OF ANALYSIS

Page 2 of 2

### Standard Curve Data:

hu IL-2 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1520	442,000	1.2
380	124,000	1.8
95.0	32,800	1.6
23.8	8,340	2.1
5.94	2,160	1.6
1.48	580	2.1
0.371	187	2.5
0.00	59	5.1


### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-2 Plate	Z0046677	2-8°C	31 Mar 2020
Anti-hu IL-2 Antibody	D0081036	2-8°C	29 Feb 2020
Proinflammatory Panel 1 (human) Calibrator Blend	A0080139	2-8°C	31 Jul 2022
Diluent 2	M0020098	≤ -10°C	28 Feb 2020
Diluent 3	M0030078	≤ -10°C	31 Jan 2020
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>	Dawn Stollar	Quality Assurance		29 NOV 2017