

**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:** K0080770  
**Expiration Date:** 30 Apr 2019

**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	1540	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.

**Standard Curve Data:**

Hu IL-2 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1540	443,000	1.4
385	126,000	4.0
96.3	33,900	2.7
24.1	8,680	1.0
6.02	2,200	1.9
1.50	620	9.7
0.376	197	2.9
0.00	59	5.1

## CERTIFICATE OF ANALYSIS

Page 2 of 2

### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-2 Plate	Z0046500	2-8°C	31 Oct 2019
Anti-hu IL-2 antibody	D0080784	2-8°C	30 Apr 2019
Proinflammatory Panel 1 (human) Calibrator Blend	A0080123	2-8°C	31 Mar 2022
Diluent 2	M0020095	≤ -10°C	30 Sep 2019
Diluent 3	M0030076	≤ -10°C	30 Sep 2019
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), 1000mL	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	<i>Dawn Stollar</i>	28 JUN 2017