

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-15 Kit  
**Kit Catalog Numbers:** K151RD-Series  
**Kit Lot Number:** K0081089  
**Expiration Date:** 31 Dec 2020

### Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Cytokine Panel 1 (human) Control 1		Cytokine Panel 1 (human) Control 2		Cytokine Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu IL-15	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu IL-15	0.774	80% - 120%	< 20%	525	80% - 120%	< 20%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IL-15	Calibrator	842	pg/mL
		Cytokine Panel 1 (human) Control 1	288	pg/mL
		Cytokine Panel 1 (human) Control 2	46.3	pg/mL
		Cytokine Panel 1 (human) Control 3	8.87	pg/mL

Controls tested may not be the same lot as shipped.

### Standard Curve Data:

hu IL-15 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
842	308,000	2.4
211	54,700	2.0
52.6	12,300	3.3
13.2	2,920	4.0
3.29	748	2.3
0.822	237	3.6
0.206	118	5.9
0.00	81	5.1

**CERTIFICATE OF ANALYSIS**


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-15 Plate	Z0046913	2-8°C	28 Feb 2021
Anti-hu IL-15 Antibody	D0081081	2-8°C	30 Sep 2021
Cytokine Panel 1 (human) Calibrator Blend	A0080144	2-8°C	31 Aug 2022
Diluent 43	M0430027	≤ -10°C	31 Dec 2020
Diluent 3	M0030084	≤ -10°C	31 Jan 2021
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), 1000 mL	Not Kit Specific
Cytokine Panel 1 (human) Control 1	Not Kit Specific
Cytokine Panel 1 (human) Control 2	Not Kit Specific
Cytokine 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		02 NOV 2018