

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® AND V-PLEX Plus Human IL-10 Kit  
**Kit Catalog Numbers:** K151QU-Series, K156QU-Series  
**Kit Lot Number:** K0081356  
**Expiration Date:** 28 Feb 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-10	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-10	0.298	80% - 120%	< 20%	233	80% - 120%	< 20%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-10	Calibrator	378	pg/mL
		Proinflammatory Panel 1 (human) Control 1	132	pg/mL
		Proinflammatory Panel 1 (human) Control 2	24.8	pg/mL
		Proinflammatory Panel 1 (human) Control 3	3.62	pg/mL

Controls tested may not be the same lot as shipped.

### Standard Curve Data:

IL-10 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
378	180,000	1.4
94.5	47,300	2.6
23.6	12,100	1.8
5.91	3,200	1.7
1.48	826	2.8
0.369	248	3.3
0.0923	104	3.7
0.00	57	6.4

## CERTIFICATE OF ANALYSIS


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-10 Plate	Z0047208	2-8°C	31 Mar 2022
Anti-hu IL-10 Antibody	D0081134	2-8°C	31 Mar 2022
Proinflammatory Panel 1 (human) Calibrator Blend	A0080186	2-8°C	30 Nov 2023
Diluent 2	M0020109	≤ -10°C	28 Feb 2022
Diluent 3	M0030092	≤ -10°C	31 Mar 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>	Carissa Hubbard	Quality		10 DEC 2019