

**CERTIFICATE OF ANALYSIS**

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-6 Kit  
**Kit Catalog Numbers:** K151QX-Series  
**Kit Lot Number:** K0081023  
**Expiration Date:** 30 Nov 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-6	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-6	0.633	80% - 120%	< 20%	488	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	hu IL-6	Calibrator	732	pg/mL
		Proinflammatory Panel 1 (human) Control 1	264	pg/mL
		Proinflammatory Panel 1 (human) Control 2	41.2	pg/mL
		Proinflammatory Panel 1 (human) Control 3	8.24	pg/mL

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

hu IL-6 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
732	617,000	14.0
183	154,000	3.2
45.8	33,700	5.1
11.4	8,070	2.5
2.86	2,060	3.5
0.715	570	2.8
0.179	192	3.3
0.00	72	5.8

## CERTIFICATE OF ANALYSIS

Page 2 of 2


### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-6 Plate	Z0046844	2-8°C	30 Nov 2020
Anti-hu IL-6 Antibody	D0081006	2-8°C	30 Nov 2019
Proinflammatory Panel 1 (human) Calibrator Blend	A0080158	2-8°C	31 May 2023
Diluent 2	M0020099	≤ -10°C	31 Mar 2020
Diluent 3	M0030082	≤ -10°C	31 Oct 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), 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		03 JUL 2018