

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-8 Kit  
**Kit Catalog Numbers:** K156RA-Series, K151RA-Series  
**Kit Lot Number:** K0081502  
**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-8	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-8	0.591	80% - 120%	< 20%	375	80% - 120%	< 20%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-8	Calibrator	560	pg/mL
		Proinflammatory Panel 1 (human) Control 1	220	pg/mL
		Proinflammatory Panel 1 (human) Control 2	33.2	pg/mL
		Proinflammatory Panel 1 (human) Control 3	5.42	pg/mL

Controls tested may not be the same lot as shipped.

### Standard Curve Data:

IL-8 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
560	300,000	2.2
140	75,400	2.2
35.0	19,000	1.6
8.75	4,850	2.1
2.19	1,280	2.0
0.547	368	2.7
0.137	141	2.8
0.00	68	6.6

## CERTIFICATE OF ANALYSIS


### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-8 Plate	Z0047347	2-8°C	30 Nov 2022
Anti-hu IL-8 Antibody	D0081142	2-8°C	28 Feb 2022
Proinflammatory Panel 1 (human) Calibrator Blend	A0080217	2-8°C	30 Jun 2024
Diluent 2	M0020113	≤ -10°C	30 Nov 2022
Diluent 3	M0030100	≤ -10°C	31 Oct 2022
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
MSD Wash Buffer (20X), 100mL	Not Kit Specific
Plate Seals	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>	Karen Hamilla	Quality		20 JUL 2020