

**CERTIFICATE OF ANALYSIS**

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-8 Kit

**Kit Catalog Numbers:** K151RA-Series, K156RA-Series

**Kit Lot Number:** K0081530

**Expiration Date:** 30 Apr 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	592	pg/ml
		Proinflammatory Panel 1(human) Control 1	233	pg/ml
		Proinflammatory Panel 1(human) Control 2	36.8	pg/ml
		Proinflammatory Panel 1(human) Control 3	6.29	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
592	375,000	3.9
148	92,000	2.4
37.0	22,600	5.0
9.25	5,650	1.2
2.31	1,420	1.8
0.578	409	2.7
0.145	157	4.4
0.00	73	4.8

## CERTIFICATE OF ANALYSIS

Page 2 of 2


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-8 Plate	Z0047373	2-8°C	31 Jan 2023
Anti-hu IL-8 Antibody	D0081168	2-8°C	30 Apr 2022
Proinflammatory Panel 1(human) Calibrator Blend	A0080229	2-8°C	30 Sep 2024
Diluent 2	M0020113	≤ -10°C	30 Nov 2022
Diluent 3	M0030101	≤ -10°C	31 Dec 2022
MSD Read Buffer T (4X)	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>	Karen Hamilla	Quality		28 OCT 2020