

**CERTIFICATE OF ANALYSIS**

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-8 Kit  
**Kit Catalog Numbers:** K151RA-Series  
**Kit Lot Number:** K0081099  
**Expiration Date:** 31 Jan 2021

**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-8	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-8	0.591	80% - 120%	< 20%	375	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	hu IL-8	Calibrator	621	pg/mL
		Proinflammatory Panel 1 (human) Control 1	235	pg/mL
		Proinflammatory Panel 1 (human) Control 2	36.2	pg/mL
		Proinflammatory Panel 1 (human) Control 3	5.93	pg/mL

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

hu IL-8 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
621	351,000	1.4
155	80,800	2.4
38.8	19,500	1.2
9.70	4,640	2.4
2.43	1,220	2.0
0.606	347	3.2
0.152	138	2.9
0.00	68	5.4

**Components:**



## CERTIFICATE OF ANALYSIS

Page 2 of 2

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-8 Plate	Z0046925	2-8°C	31 Mar 2021
Anti-hu IL-8 Antibody	D0081097	2-8°C	30 Nov 2021
Proinflammatory Panel 1 (human) Calibrator Blend	A0080178	2-8°C	31 Aug 2023
Diluent 2	M0020102	≤ -10°C	28 Feb 2021
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), 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>	Dawn Stollar	Quality Assurance		22 OCT 2018