

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: K0081015
Expiration Date: 31 Dec 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-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	561	pg/mL
		Proinflammatory Panel 1 (human) Control 1	233	pg/mL
		Proinflammatory Panel 1 (human) Control 2	36.9	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
561	276,000	2.4
140	68,600	1.3
35.1	17,100	3.1
8.77	4,280	1.6
2.19	1,150	1.3
0.548	351	2.0
0.137	153	3.2
0.00	95	28.0

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-8 Plate	Z0046832	2-8°C	31 Oct 2020
Anti-hu IL-8 Antibody	D0081012	2-8°C	31 Dec 2019
Proinflammatory Panel 1 (human) Calibrator Blend	A0080158	2-8°C	31 May 2023
Diluent 2	M0020099	≤ -10°C	31 Mar 2020
Diluent 3	M0030081	≤ -10°C	30 Sep 2020
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

CERTIFICATE OF ANALYSIS

Page 2 of 2

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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	25 JUN 2018