

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-17A Kit
Kit Catalog Numbers: K151RF-Series; K156RF-Series
Kit Lot Number: K0081295
Expiration Date: 31 Dec 2021

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Cytokine Panel 1 (human) Control 1		Cytokine Panel 1 (human) Control 2		Cytokine Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-17A	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-17A	3.19	80% - 120%	< 20%	3650	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-17A	Calibrator	5770	pg/mL
		Cytokine Panel 1 (human) Control 1	2590	pg/mL
		Cytokine Panel 1 (human) Control 2	374	pg/mL
		Cytokine Panel 1 (human) Control 3	58.6	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IL-17A - Spot 1		
Conc. pg/mL	ECL Signal	%CV
5770	558,000	2.3
1440	144,000	3.2
361	35,500	5.1
90.2	8,840	3.8
22.5	2,240	3.3
5.63	613	5.7
1.41	189	3.6
0.00	56	9.2

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-17A Plate	Z0047149	2-8°C	31 Jan 2022
Anti-hu IL-17A Antibody	D0081187	2-8°C	30 Apr 2022
Cytokine Panel 1 (human) Calibrator Blend	A0080227	2-8°C	31 Oct 2024
Diluent 43	M0430033	≤ -10°C	31 Jan 2022
Diluent 3	M0030090	≤ -10°C	31 Dec 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
Cytokine Panel 1 (human) Control 1	Not Kit Specific
Cytokine Panel 1 (human) Control 2	Not Kit Specific
Cytokine 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	Karen Hamilla	Quality		29 JAN 2020