

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-17A/F Kit
Kit Catalog Numbers: K151WN-Series
Kit Lot Number: K0081406
Expiration Date: 31 Jan 2022

Acceptance Data/ Percent Recovery Specifications:

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

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-17A/F	7.57	80% - 120%	< 20%	3900	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-17A/F	Calibrator	6140	pg/mL
		Cytokine Panel 2 (human) Control 1	5550	pg/mL
		Cytokine Panel 2 (human) Control 2	636	pg/mL
		Cytokine Panel 2 (human) Control 3	84.2	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IL-17A/F - Spot 1		
Conc. pg/mL	ECL Signal	%CV
6140	401,000	3.2
1540	94,100	1.5
384	23,300	1.9
95.9	5,920	3.0
24.0	1,560	1.9
6.00	473	2.1
1.50	190	2.9
0.00	98	3.6

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-17A/F Plate	Z0047253	2-8°C	30 Jun 2022
Anti-hu IL-17A/F Antibody	D0081425	2-8°C	30 Jun 2023
Cytokine Panel 2 (human) Calibrator Blend	A0080116	2-8°C	31 Jan 2022
Diluent 43	M0430036	≤ -10°C	31 May 2022
Diluent 3	M0030094	≤ -10°C	31 May 2022
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 2 (human) Control 1	Not Kit Specific
Cytokine Panel 2 (human) Control 2	Not Kit Specific
Cytokine Panel 2 (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		10 MAR 2020