

**CERTIFICATE OF ANALYSIS**

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-17A Gen. B Kit  
**Kit Catalog Numbers:** K151WM-Series  
**Kit Lot Number:** K0081517  
**Expiration Date:** 31 Mar 2022

**Acceptance Data/ Percent Recovery Specifications:**

Spot	Name	Calibrator		TH17 Panel 1 (human) Control 1		TH17 Panel 1 (human) Control 2		TH17 Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-17A	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	5.86	80% - 120%	< 20%	1950	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	IL-17A	Calibrator	2570	pg/mL
		TH17 Panel 1 (human) Control 1	2860	pg/mL
		TH17 Panel 1 (human) Control 2	403	pg/mL
		TH17 Panel 1 (human) Control 3	57.1	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
2570	158,000	5.0
643	37,500	3.8
161	9,190	3.9
40.2	2,340	3.5
10.0	658	3.6
2.51	219	3.6
0.627	110	5.8
0.00	75	5.4

## CERTIFICATE OF ANALYSIS

Page 2 of 2


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-17A Gen. B Plate	Z0047354	2-8°C	31 Dec 2022
Anti-hu IL-17A Antibody	D0081135	2-8°C	30 Apr 2022
TH17 Panel 1 (human) Calibrator Blend	A0080232	2-8°C	31 Oct 2024
Diluent 43	M0430035	≤ -10°C	31 Mar 2022
Diluent 3	M0030092	≤ -10°C	31 Mar 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), 100 mL	Not Kit Specific
TH17 Panel 1 (human) Control 1	Not Kit Specific
TH17 Panel 1 (human) Control 2	Not Kit Specific
TH17 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		17 SEP 2020