



## RFE MEMO

Page 1 of 1

**Product Description(s):** V-PLEX® and V-PLEX Plus Human IL-17A Gen. B Kit


**Catalog Number(s) / Size(s):** K151WM-Series

**Lot Number:** K0081490

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0081490 been extended to 30 APR 2023.

The expiration date for the following kit component has been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Human IL-17A Gen. B Plate	Z0047331	31 OCT 2022	30 APR 2023
Anti-hu IL-17A Antibody	D0081135	30 APR 2022	30 APR 2023
Diluent 43	M0430036	31 MAY 2022	30 APR 2023
Diluent 3	M0030096	30 JUN 2022	30 APR 2023

Function	Name	Signature	Date
Quality	Karen Hamilla		11 APR 2022

This memo supersedes all other statements regarding date of expiration for this product.

## 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:** K0081490  
**Expiration Date:** 30 Apr 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 Gen. B	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 Gen. B	5.86	80% - 120%	< 20%	1950	80% - 120%	< 20%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-17A Gen. B	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 Gen. B - Spot 1		
Conc. pg/mL	ECL Signal	%CV
2570	102,000	2.9
643	22,900	3.7
161	5,690	2.9
40.2	1,370	4.1
10.0	430	3.3
2.51	177	2.5
0.627	115	2.3
0.00	92	3.5

**CERTIFICATE OF ANALYSIS**

**Components:**


Description	Lot Number	Storage Temperature	Expiration Date
Human IL-17 A Gen. B Plate	Z0047331	2-8°C	31 Oct 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	M0430036	≤ -10°C	31 May 2022
Diluent 3	M0030096	≤ -10°C	30 Jun 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
<b>Review/Approval</b>	Karen Hamilla	Quality		19 JUN 2020