



## RFE MEMO

Page 1 of 1

**Product Description(s):** V-PLEX® and V-PLEX Plus Human IL-22 Kit


**Catalog Number(s) / Size(s):** K151WV - Series

**Lot Number:** K0081412

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0081412 has been extended to 28 FEB 2023.

The expiration dates for the following kit components have been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Human IL-22 Plate	Z0047256	31 JUL 2022	28 FEB 2023
Anti-hu IL-22 Antibody	D0081137	28 FEB 2022	28 FEB 2023
Diluent 43	M0430036	31 MAY 2022	28 FEB 2023
Diluent 3	M0030094	31 MAY 2022	28 FEB 2023

Function	Name	Signature	Date
Quality	Karen Hamilla		21 JAN 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-22 Kit

**Kit Catalog Numbers:** K151WV-Series

**Kit Lot Number:** K0081412

**Expiration Date:** 28 Feb 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-22	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-22	2.78	80% - 120%	< 20%	325	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	IL-22	Calibrator	527	pg/mL
		TH17 Panel 1 (human) Control 1	525	pg/mL
		TH17 Panel 1 (human) Control 2	84.7	pg/mL
		TH17 Panel 1 (human) Control 3	17.6	pg/mL

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

IL-22 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
527	112,000	2.5
132	28,700	1.4
32.9	7,110	2.1
8.23	1,910	2.6
2.06	563	2.4
0.515	236	1.8
0.129	153	1.9
0.00	123	4.1

## CERTIFICATE OF ANALYSIS

Page 2 of 2


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-22 Plate	Z0047256	2-8°C	31 Jul 2022
Anti-hu IL-22 Antibody	D0081137	2-8°C	28 Feb 2022
TH17 Panel 1 (human) Calibrator Blend	A0080232	2-8°C	31 Oct 2024
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
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		11 MAR 2020