

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-22 Kit  
**Kit Lot Number:** K0082020  
**Expiration Date:** 31 Dec 2023  
**Kit Catalog Numbers:** K151WV-Series

**Precision and Accuracy 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 Calibrator:**


Spot	Name	Conc.	Units
1	IL-22	504	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-22 Plate	Z0047696	2-8°C	31 Oct 2024
Anti-hu IL-22 Antibody	D0081522	2-8°C	31 Dec 2023
TH17 Panel 1 (human) Calibrator Blend	A0080303	2-8°C	31 Aug 2026
Diluent 43	M0430047	≤ -10°C	31 May 2024
Diluent 3	M0030114	≤ -10°C	30 Sep 2024

Additional components may be provided to execute the assay. Additional components are not lot-specific.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		14 JUL 2022