

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX® and V-PLEX Plus Human IL-17A Gen. B Kit
Lot Number: K0082451
Expiration Date: 30 NOV 2025
Catalog Number: K151WM-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IL-17A Antibody	D0081862	2-8°C	30 NOV 2025
Human IL-17A Gen. B SECTOR Plate	Z0047976	2-8°C	30 SEP 2026
TH17 Panel 1 (human) Calibrator Blend	A0080366	2-8°C	31 MAY 2028
Diluent 3	M0030123	≤ -10°C	31 JAN 2026
Diluent 43	M0430053	≤ -10°C	31 JAN 2026

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

Precision, Accuracy, and Sensitivity 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 Calibrator:

Spot	Name	Conc.	Units
1	IL-17A Gen.B	3230	pg/mL

Dilution to top of Curve: 1X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		02 MAY 2024