



EXTENSION 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: K0081936

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0081936 has been extended to 28 FEB 2025.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Human IL-17A Gen. B SECTOR Plate	Z0047642	30 JUN 2024	28 FEB 2025
Anti-hu IL-17A Antibody	D0081613	31 JUL 2024	28 FEB 2025
Diluent 43	M0430046	29 FEB 2024	28 FEB 2025
Diluent 3	M0030110	31 MAR 2024	28 FEB 2025

Function	Name	Signature	Date
Quality	Karen Hamilla		08 DEC 2023

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

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-Plex® and V-Plex Plus Human IL-17A Gen. B Kit
Kit Lot Number: K0081936
Expiration Date: 29 Feb 2024
Kit Catalog Numbers: K151WM-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-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	3140	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-17A Gen. B SECTOR Plate	Z0047642	2-8°C	30 Jun 2024
Anti-hu IL-17A Antibody	D0081613	2-8°C	31 Jul 2024
TH17 Panel 1 (human) Calibrator Blend	A0080241	2-8°C	28 Feb 2025
Diluent 43	M0430046	≤ -10°C	29 Feb 2024
Diluent 3	M0030110	≤ -10°C	31 Mar 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		03 MAY 2022