



Extension Memo

Page 1 of 1

Product Description: V-PLEX® and V-PLEX PLUS Human TARC Gen. B Kit
Catalog Number(s): K151ANK-Series, K156ANK-Series
Lot Number: K0082352
New Expiration Date: 31 Jan 2026

Based on additional testing, the original date of expiration for this kit has been extended by 6 months.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Diluent 57	M0570011	31 Jul 2025	31 Jan 2026

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		08 Aug 2025



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX PLUS Human TARC Gen. B Kit
Lot Number: K0082352
Expiration Date: 31 JUL 2025
Catalog Number: K151ANK-Series; K156ANK-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Chemokine Panel 1 (human) Gen. B Calibrator Blend	A0080377	2-8°C	31 AUG 2028
Human TARC Antibody	D0082132	2-8°C	31 JAN 2027
Human TARC Gen. B Plate	Z0047903	2-8°C	28 FEB 2026
Diluent 3	M0030123	≤ -10°C	31 JAN 2026
Diluent 57	M0570011	≤ -10°C	31 JUL 2025

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

Precision, Accuracy, and Sensitivity Specifications:

Spot	Name	Calibrator		Chemokine Panel 1 (human) Gen. B Control 1		Chemokine Panel 1 (human) Gen. B Control 2		Chemokine Panel 1 (human) Gen. B Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	TARC/CCL17	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	TARC/CCL17	2.14	80% - 120%	< 20%	481	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	TARC/CCL17	744	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		30 MAY 2024