



EXTENSION MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX Plus Human TARC Kit


Catalog Number(s) / Size(s): K151NT-Series; K156NT-Series

Lot Number: K0082371

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0082371 has been extended to 30 NOV 2024 .

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

Description	Lot Number	Original Expiration Date	New Expirati
Human TARC Plate	Z0047539	30 NOV 2023	30 NOV 2024
Diluent 43	M0430049	31 OCT 2024	30 NOV 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		09 OCT 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 TARC Kit
Kit Lot Number: K0082371
Expiration Date: 30 Nov 2023
Kit Catalog Numbers: K151NT-Series; K156NT-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human TARC Plate	Z0047539	2-8°C	30 Nov 2023
Anti-hu TARC Antibody	D0081681	2-8°C	31 Jul 2025
Chemokine Panel 1 (human) Calibrator Blend	A0080306	2-8°C	30 Sep 2026
Diluent 3	M0030120	≤ -10°C	31 Jul 2025
Diluent 43	M0430049	≤ -10°C	31 Oct 2024

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

Precision and Accuracy Specifications:

		Calibrator		Chemokine Panel 1 (human) Control 1		Chemokine Panel 1 (human) Control 2		Chemokine Panel 1 (human) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	TARC	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%


		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	TARC	3.32	80% - 120%	< 20%	1120	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	TARC	2080	pg/mL

Dilution to Top of Curve = 1X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		09 OCT 2023