



## CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX® and V-PLEX Plus Human TARC Kit  
**Lot Number:** K0082571  
**Expiration Date:** 30 JUN 2026  
**Catalog Number:** K151NT-Series, K156NT-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu TARC Antibody	D0081968	2-8°C	30 JUN 2026
Chemokine Panel 1 (human) Calibrator Blend	A0080388	2-8°C	31 MAR 2029
Human TARC SECTOR Plate	Z0048032	2-8°C	31 MAR 2027
Diluent 3	M0030128	≤ -10°C	31 OCT 2026
Diluent 43	M0430056	≤ -10°C	30 SEP 2026

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

Spot	Name	LLOQ			ULOQ		
		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	1750	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
<b>Review/Approval</b>	Karen Hamilla	Quality		19 DEC 2024