



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Human TARC Kit
Lot Number: K0082711
Expiration Date: 30 SEP 2026
Catalog Number: K151NT-Series; K156NT-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu TARC Antibody	D0082174	2-8°C	30 JUN 2027
Chemokine Panel 1 (human) Calibrator Blend	A0080388	2-8°C	31 MAR 2029
Human TARC 96-Well Small Spot 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
Review/Approval	Karen Hamilla	Quality		01 MAY 2025