



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Human TARC Kit**Lot Number:** K0082777**Expiration Date:** 28 FEB 2027**Catalog Number:** K151NT-Series, K156NT-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu TARC Antibody	D0082475	2-8°C	30 SEP 2028
Chemokine Panel 1 (human) Calibrator Blend	A0080432	2-8°C	30 APR 2030
Human TARC SECTOR Plate	Z0048165	2-8°C	29 FEB 2028
Diluent 3	M0030141	≤ -10°C	30 NOV 2027
Diluent 43	M0430058	≤ -10°C	28 FEB 2027

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	1780	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		17 OCT 2025