

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Human TARC Kit  
**Kit Lot Number:** K0082383  
**Expiration Date:** 30 Sep 2025  
**Kit Catalog Numbers:** K151NT-Series; K156NT-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human TARC Plate	Z0047919	2-8°C	31 Mar 2026
Anti-hu TARC Antibody	D0081767	2-8°C	30 Sep 2025
Chemokine Panel 1 (human) Calibrator Blend	A0080364	2-8°C	31 May 2028
Diluent 3	M0030123	≤ -10°C	31 Jan 2026
Diluent 43	M0430053	≤ -10°C	31 Jan 2026

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

**Precision and Accuracy 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%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 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	1690	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		13 Nov 2023