



Meso Scale Diagnostics, LLC.

### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human TARC Kit  
**Kit Lot Number:** K0082328  
**Expiration Date:** 31 Jul 2025  
**Kit Catalog Numbers:** K151NT-Series; K156NT-Series

#### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human TARC Plate	Z0047889	2-8°C	31 Dec 2025
Anti-hu TARC Antibody	D0081681	2-8°C	31 Jul 2025
Chemokine Panel 1 (human) Calibrator Blend	A0080364	2-8°C	31 May 2028
Diluent 3	M0030123	≤ -10°C	31 Jan 2026
Diluent 43	M0430052	≤ -10°C	31 Dec 2025

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		14 Sep 2023