

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human TARC Kit  
**Kit Lot Number:** K0081592  
**Expiration Date:** 30 Jun 2022  
**Kit Catalog Numbers:** K151NT-Series, K156NT-Series

### Precision and Accuracy Specifications:

		Calibrator		Chemokine Panel 1 (human) Control 1		Chemokine Panel 1 (human) Control 2		Chemokine Panel 1 (human) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	TARC	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

		LLOQ			ULOQ		
Spot	Name	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	1845	pg/mL

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human TARC plate	Z0047377	2-8°C	28 Feb 2023
Anti-hu TARC Antibody	D0081261	2-8°C	31 Aug 2022
Chemokine Panel 1 (human) Calibrator Blend	A0080266	2-8°C	31 Aug 2025
Diluent 43	M0430037	≤ -10°C	31 Aug 2022
Diluent 3	M0030096	≤ -10°C	30 Jun 2022

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

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		19 NOV 2020