

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human TARC Kit
Kit Lot Number: K0081626
Expiration Date: 31 Aug 2022
Kit Catalog Numbers: K151NT-Series, K156NT-Series

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	1845	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human TARC Plate	Z0047444	2-8°C	30 Jun 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	M0430043	≤ -10°C	31 Jul 2023
Diluent 3	M0030104	≤ -10°C	30 Jun 2023

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	Xiomara Talavera	Quality		19 APR 2021