

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human TARC Kit

Kit Catalog Numbers: K151NT-Series, K156NT-Series

Kit Lot Number: K0081428

Expiration Date: 31 Oct 2021

Acceptance Data/ Percent Recovery 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 Material Tested :

Spot	Name	Item	Conc	Units
1	TARC	Calibrator	1870	pg/mL
		Chemokine Panel 1 (human) Control 1	1300	pg/mL
		Chemokine Panel 1 (human) Control 2	216	pg/mL
		Chemokine Panel 1 (human) Control 3	37	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

TARC - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1870	301,000	5.2
468	68,500	5.6
117	16,700	5.0
29.2	4,300	6.8
7.30	1,150	4.5
1.83	331	3.3
0.457	125	4.3
0.00	55	7.1

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human TARC plate	Z0047265	2-8°C	31 Jul 2022
Anti-hu TARC Antibody	D0081090	2-8°C	31 Oct 2021
Chemokine Panel 1 (human) Calibrator Blend	A0080204	2-8°C	30 Apr 2024
Diluent 43	M0430036	≤ -10°C	31 May 2022
Diluent 3	M0030096	≤ -10°C	30 Jun 2022
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Plate Seals	Not Kit Specific
Wash Buffer (20X), 100mL	Not Kit Specific
Chemokine Panel 1 (human) Control 1	Not Kit Specific
Chemokine Panel 1 (human) Control 2	Not Kit Specific
Chemokine Panel 1 (human) Control 3	Not Kit Specific

Statement:

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		24 MAR 2020