

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human TARC Kit
Kit Catalog Numbers: K151NT-Series
Kit Lot Number: K0080980
Expiration Date: 31 Jul 2019

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	hu TARC	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu TARC	3.32	80% - 120%	< 20%	1120	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu TARC	Calibrator	1600	pg/mL
		Chemokine Panel 1 (human) Control 1	1390	pg/mL
		Chemokine Panel 1 (human) Control 2	210	pg/mL
		Chemokine Panel 1 (human) Control 3	36.5	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu TARC - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1600	388,000	2.9
400	95,900	2.5
100	23,600	3.5
25.0	5,920	2.6
6.25	1,610	4.2
1.56	446	2.2
0.391	165	3.1
0.00	73	4.5

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human TARC Plate	Z0046790	2-8°C	31 Aug 2020
Anti-hu TARC Antibody	D0080900	2-8°C	31 Jul 2019
Chemokine Panel 1 (human) Calibrator Blend	A0080159	2-8°C	31 Dec 2022
Diluent 43	M0430021	≤ -10°C	30 Sep 2019
Diluent 3	M0030076	≤ -10°C	30 Sep 2019
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Wash Buffer (20X), 100mL	Not Kit Specific
Plate Seals	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	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	04 JUN 2018