

CERTIFICATE OF ANALYSIS

Page 1 of 2

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

Kit Catalog Numbers: K156NT-Series

Kit Lot Number: K0080853

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	NHP TARC	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

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

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	NHP TARC	Calibrator	1890	pg/mL
		Chemokine Panel 1 (human) Control 1	1130	pg/mL
		Chemokine Panel 1 (human) Control 2	197	pg/mL
		Chemokine Panel 1 (human) Control 3	34.2	pg/mL

Controls tested may not be the same lot as shipped.

The calibrator and controls for this assay contain a human recombinant cytokine calibrator sequence that is highly homologous to the non-human primate cytokine. This assay has been verified to cross-react with homologs of rhesus and cynomolgus monkeys. See Product Insert for experimental data.

Standard Curve Data:

NHP TARC - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1890	391,000	5.2
473	111,000	4.6
118	28,400	4.1
29.5	7,050	4.6
7.38	1,980	4.4
1.85	554	5.1
0.461	187	3.5
0.00	60	5.6

Components:

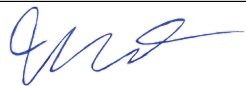
CERTIFICATE OF ANALYSIS

Description	Lot Number	Storage Temperature	Expiration Date
Human TARC Plate	Z0046583	2-8°C	31 Dec 2019
Anti-hu TARC Antibody	D0080900	2-8°C	31 Jul 2019
Chemokine Panel 1 (human) Calibrator Blend	A0080122	2-8°C	28 Feb 2022
Diluent 43	M0430023	≤ -10°C	31 Dec 2019
Diluent 3	M0030077	≤ -10°C	30 Nov 2019
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	Jessica Drolet	Quality Assurance		03 May 2018