

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: K0081178
Expiration Date: 28 Feb 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	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	1800	pg/mL
		Chemokine Panel 1 (human) Control 1	1350	pg/mL
		Chemokine Panel 1 (human) Control 2	213	pg/mL
		Chemokine Panel 1 (human) Control 3	36	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
1800	345,000	4.3
450	89,200	5.7
113	22,400	5.7
28.1	5,700	4.9
7.03	1,460	6.9
1.76	414	5.5
0.439	151	5.6
0.00	67	5.9

CERTIFICATE OF ANALYSIS

Components:

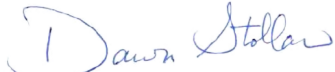
Description	Lot Number	Storage Temperature	Expiration Date
Human TARC Plate	Z0047008	2-8°C	30 Jun 2021
Anti-hu TARC Antibody	D0081163	2-8°C	28 Feb 2022
Chemokine Panel 1 (human) Calibrator Blend	A0080183	2-8°C	31 Oct 2023
Diluent 43	M0430029	≤ -10°C	30 Apr 2021
Diluent 3	M0030085	≤ -10°C	28 Feb 2021
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), 100 mL	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		18 MAR 2019