

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus NHP IL-7 Kit

Kit Catalog Numbers: K156RC-Series

Kit Lot Number: K0080971

Expiration Date: 31 Jul 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Cytokine Panel 1(human) Control 1		Cytokine Panel 1(human) Control 2		Cytokine Panel 1(human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	NHP IL-7	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	NHP IL-7	1.37	80% - 120%	< 20%	563	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	NHP IL-7	Calibrator	1010	pg/mL
		Cytokine Panel 1(human) Control 1	330	pg/mL
		Cytokine Panel 1(human) Control 2	60	pg/mL
		Cytokine Panel 1(human) Control 3	10	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 IL-7 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1010	241,000	2.1
253	58,200	2.1
63.1	14,500	1.9
15.8	3,710	2.6
3.95	1,010	1.4
0.986	309	2.1
0.00	133	2.6

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-7 Plate	Z0046777	2-8°C	31 Aug 2020
Anti-hu IL-7 Antibody	D0080888	2-8°C	31 Jul 2019
Cytokine Panel 1 (human) Calibrator Blend	A0080108	2-8°C	30 Nov 2021
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
Cytokine Panel 1 (human) Control 1	Not Kit Specific
Cytokine Panel 1 (human) Control 2	Not Kit Specific
Cytokine 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>	10 APR 2018