

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: K0081184
Expiration Date: 28 Feb 2021

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	0.546	80% - 120%	< 20%	563	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	NHP IL-7	Calibrator	846	pg/mL
		Cytokine Panel 1 (human) Control 1	346	pg/mL
		Cytokine Panel 1 (human) Control 2	58	pg/mL
		Cytokine Panel 1 (human) Control 3	7.79	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
846	193,000	2.5
212	46,100	2.1
52.9	11,300	2.8
13.2	2,770	3.1
3.30	729	2.6
0.826	238	2.7
0.207	123	4.9
0.00	84	6.0

CERTIFICATE OF ANALYSIS

Page 2 of 2

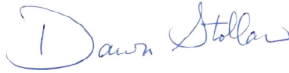
Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-7 Plate	Z0047016	2-8°C	30 Jun 2021
Anti-hu IL-7 Antibody	D0081079	2-8°C	30 Sep 2021
Cytokine Panel 1 (human) Calibrator Blend	A0080170	2-8°C	30 Jun 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), 100mL	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		05 MAR 2019