

## 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:** K0081163  
**Expiration Date:** 30 Apr 2020

### 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	819	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
819	193,000	2.6
205	47,400	1.9
51.2	11,900	1.8
12.8	2,970	1.2
3.20	814	1.3
0.800	250	1.9
0.200	114	4.3
0.00	70	4.0

**CERTIFICATE OF ANALYSIS**


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-7 Plate	Z0046990	2-8°C	31 May 2021
Anti-hu IL-7 Antibody	D0081055	2-8°C	30 Apr 2020
Cytokine Panel 1(human) Calibrator Blend	A0080063	2-8°C	31 Jul 2020
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	
MSD Wash Buffer (20X), 1000mL	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
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		14 JAN 2019