

**CERTIFICATE OF ANALYSIS**

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus NHP IL-15 Kit

**Kit Catalog Numbers:** K156RD-Series

**Kit Lot Number:** K0080941

**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-15	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-15	1.40	80% - 120%	< 20%	525	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	NHP IL-15	Calibrator	842	pg/mL
		Cytokine Panel 1 (human) Control 1	241	pg/mL
		Cytokine Panel 1 (human) Control 2	36.5	pg/mL
		Cytokine Panel 1 (human) Control 3	8	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-15 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
842	272,000	2.1
211	47,800	3.8
52.6	11,100	2.5
13.2	2,860	3.1
3.29	748	2.0
0.822	237	2.8
0.206	107	3.7
0.00	64	5.9

**CERTIFICATE OF ANALYSIS**

**Components:**


Description	Lot Number	Storage Temperature	Expiration Date
Human IL-15 Plate	Z0046720	2-8°C	30 Jun 2020
Anti-hu IL-15 Antibody	D0080890	2-8°C	31 Jul 2019
Cytokine Panel 1 (human) Calibrator Blend	A0080144	2-8°C	31 Aug 2022
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**

Plate Seals	Not Kit Specific
Wash Buffer (20x), 1000mL	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		31 JAN 2018