

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: K0080988

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	757	pg/mL
		Cytokine Panel 1 (human) Control 1	280	pg/mL
		Cytokine Panel 1 (human) Control 2	50	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-15 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
757	279,000	1.7
189	48,900	2.0
47.3	11,300	1.8
11.8	2,790	1.8
2.96	763	5.4
0.739	236	3.2
0.185	105	4.8
0.00	63	5.2

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:


Description	Lot Number	Storage Temperature	Expiration Date
Human IL-15 Plate	Z0046798	2-8°C	30 Sep 2020
Anti-hu IL-15 Antibody	D0080890	2-8°C	31 Jul 2019
Cytokine Panel 1(Human) Calibrator Blend	A0080063	2-8°C	31 Jul 2020
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
Review/Approval	Dawn Stollar	Quality Assurance		24 MAY 2018