

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: K0081135
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-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	0.774	80% - 120%	< 20%	525	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	NHP IL-15	Calibrator	789	pg/mL
		Cytokine Panel 1 (human) Control 1	288	pg/mL
		Cytokine Panel 1 (human) Control 2	46.3	pg/mL
		Cytokine Panel 1 (human) Control 3	8.87	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.

CERTIFICATE OF ANALYSIS
Standard Curve Data:

NHP IL-15 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
789	294,000	2.8
197	56,300	2.0
49.3	12,100	2.4
12.3	3,020	4.2
3.08	777	3.7
0.771	237	3.1
0.193	107	3.4
0.00	68	6.7


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-15 Plate	Z0046964	2-8°C	31 Mar 2021
Anti-hu IL-15 Antibody	D0081081	2-8°C	30 Sep 2021
Cytokine Panel 1 (human) Calibrator Blend	A0080121	2-8°C	28 Feb 2022
Diluent 43	M0430028	≤ -10°C	31 Jan 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		13 NOV 2018