

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus NHP IL-12/IL-23p40 Kit

Kit Catalog Numbers: K156RJ-Series

Kit Lot Number: K0081179

Expiration Date: 31 Mar 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-12/IL-23p40	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	NHP IL-12/IL-23p40	1.32	80% - 120%	< 20%	2250	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	NHP IL-12/IL-23p40	Calibrator	3650	pg/mL
		Cytokine Panel 1 (human) Control 1	1430	pg/mL
		Cytokine Panel 1 (human) Control 2	204	pg/mL
		Cytokine Panel 1 (human) Control 3	32.6	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-12/IL-23p40 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
3650	301,000	3.8
913	80,200	2.9
228	20,500	2.6
57.0	5,140	2.9
14.3	1,370	2.7
3.56	446	3.5
0.891	212	5.0
0.00	143	4.9

CERTIFICATE OF ANALYSIS

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-12/IL-23p40 Plate	Z0047009	2-8°C	30 Jun 2021
Anti-hu IL-12/23p40 Antibody	D0081040	2-8°C	31 Mar 2020
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	<i>Dawn Stollar</i>	11 FEB 2019