

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: K0081084
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	1380	pg/mL
		Cytokine Panel 1 (human) Control 2	230	pg/mL
		Cytokine Panel 1 (human) Control 3	34.7	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	276,000	1.9
913	71,000	2.2
228	17,800	1.4
57.0	4,480	1.2
14.3	1,200	1.4
3.56	405	2.9
0.891	210	5.1
0.00	145	4.6

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-12/IL-23p40 Plate	Z0046908	2-8°C	28 Feb 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	M0430027	≤ -10°C	31 Dec 2020
Diluent 3	M0030084	≤ -10°C	31 Jan 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		08 OCT 2018