

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-22 Kit
Kit Catalog Numbers: K151WV-Series
Kit Lot Number: K0081294
Expiration Date: 30 Jun 2021

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		TH17 Panel 1 (human) Control 1		TH17 Panel 1 (human) Control 2		TH17 Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-22	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-22	2.78	80% - 120%	< 20%	325	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-22	Calibrator	527	pg/mL
		TH17 Panel 1 (human) Control 1	525	pg/mL
		TH17 Panel 1 (human) Control 2	84.7	pg/mL
		TH17 Panel 1 (human) Control 3	17.6	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IL-22 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
527	121,000	1.5
132	31,100	1.8
32.9	7,870	2.1
8.23	2,020	2.2
2.06	579	2.1
0.515	221	3.3
0.129	135	3.9
0.00	105	3.9

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-22 Plate	Z0047147	2-8°C	31 Jan 2022
Anti-hu IL-22 Antibody	D0081137	2-8°C	28 Feb 2022
TH17 Panel 1 (human) Calibrator Blend	A0080232	2-8°C	31 Oct 2024
Diluent 43	M0430030	≤ -10°C	30 Jun 2021
Diluent 3	M0030087	≤ -10°C	30 Jun 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), 100 mL	Not Kit Specific
TH17 Panel 1 (human) Control 1	Not Kit Specific
TH17 Panel 1 (human) Control 2	Not Kit Specific
TH17 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	Karen Hamilla	Quality		26 NOV 2019