

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: K0081377
Expiration Date: 28 Feb 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	115,000	2.6
132	28,500	4.9
32.9	6,990	4.1
8.23	1,840	3.1
2.06	594	2.5
0.515	277	4.4
0.129	202	3.8
0.00	176	4.8

CERTIFICATE OF ANALYSIS

Components:


Description	Lot Number	Storage Temperature	Expiration Date
Human IL-22 Plate	Z0047225	2-8°C	30 Apr 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	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 Seal	Not Kit Specific
Wash Buffer (20X), 100mL	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		11 DEC 2019