

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: K0080616
Expiration Date: 31 Jan 2019

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	Hu 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	Hu IL-22	2.78	80% - 120%	< 20%	325	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Hu IL-22	Calibrator	560	pg/mL
		TH17 Panel 1 (human) Control 1	279	pg/mL
		TH17 Panel 1 (human) Control 2	47.7	pg/mL
		TH17 Panel 1 (human) Control 3	9.33	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

Hu IL-22 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
560	161,000	1.5
140	42,400	1.8
35.0	10,600	1.6
8.75	2,860	1.1
2.19	934	1.5
0.547	466	2.9
0.137	348	2.9
0.00	304	2.1



CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-22 Plate	Z0046175	2-8°C	31 Jan 2019
Anti-hu IL-22 Antibody	D0080758	2-8°C	28 Feb 2019
TH17 Panel 1 (human) Calibrator Blend	A0080101	2-8°C	30 Sep 2021
Diluent 43	M0430018	≤ -10°C	31 Jan 2019
Diluent 3	M0030073	≤ -10°C	31 Mar 2019
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	Dawn Stollar	Quality Assurance		15 FEB 2017