

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-21 Kit
Kit Catalog Numbers: K151WU-Series
Kit Lot Number: K0080615
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-21	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Hu IL-21	1.65	80% - 120%	< 20%	650	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Hu IL-21	Calibrator	1040	pg/mL
		TH17 Panel 1 (human) Control 1	590	pg/mL
		TH17 Panel 1 (human) Control 2	82.9	pg/mL
		TH17 Panel 1 (human) Control 3	14.5	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

Hu IL-21 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1040	181,000	2.9
260	50,900	1.7
65.0	12,600	2.7
16.3	3,270	1.7
4.06	860	1.6
1.02	275	3.2
0.254	114	4.0
0.00	64	5.1



CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-21 Plate	Z0046174	2-8°C	31 Jan 2019
Anti-hu IL-21 Antibody	D0080757	2-8°C	31 Jan 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), 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	Dawn Stollar	Quality Assurance		15 FEB 2017