

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-17D Kit
Kit Catalog Numbers: K151WQ - Series
Kit Lot Number: K0080673
Expiration Date: 30 Apr 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		V-PLEX Cytokine Panel 2 (human) Control 1		V-PLEX Cytokine Panel 2 (human) Control 2		V-PLEX Cytokine Panel 2 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu IL-17D	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-17D	11.2	75% - 125%	< 20%	5200	75% - 125%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IL-17D	Calibrator	7890	pg/mL
		Cytokine Panel 2 (human) Control 1	3350	pg/mL
		Cytokine Panel 2 (human) Control 2	250	pg/mL
		Cytokine Panel 2 (human) Control 3	31.5	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu IL-17D - Spot 1		
Conc. pg/mL	ECL Signal	%CV
7890	186,000	2.9
1970	51,000	3.5
493	12,900	2.2
123	3,400	1.9
30.8	918	2.7
7.71	295	2.2
1.93	129	2.5
0.00	78	3.4

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-17D Plate	Z0046230	2-8°C	31 May 2019
Anti-hu IL-17D Antibody	D0080769	2-8°C	30 Apr 2019
Cytokine Panel 2 (human) Calibrator Blend	A0080114	2-8°C	31 Dec 2021
Diluent 43	M0430020	≤ -10°C	30 Jun 2019
Diluent 3	M0030074	≤ -10°C	30 Jun 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
Cytokine Panel 2 (human) Control 1	Not Kit Specific
Cytokine Panel 2 (human) Control 2	Not Kit Specific
Cytokine Panel 2 (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	<i>Dawn Stollar</i>	21 JUL 2017