

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: K0080753
Expiration Date: 30 Apr 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Cytokine Panel 2 (human) Control 1		Cytokine Panel 2 (human) Control 2		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	7240	pg/mL
		Cytokine Panel 2 (human) Control 1	5479	pg/mL
		Cytokine Panel 2 (human) Control 2	559	pg/mL
		Cytokine Panel 2 (human) Control 3	146	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
7240	167,000	4.8
1810	44,900	3.9
453	11,000	3.9
113	2,820	4.4
28.3	781	2.7
7.07	263	2.8
1.77	128	2.3
0.00	85	6.0

CERTIFICATE OF ANALYSIS

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	A0080116	2-8°C	31 Jan 2022
Diluent 43	M0430021	≤ -10°C	30 Sep 2019
Diluent 3	M0030076	≤ -10°C	30 Sep 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)	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>	05 APR 2018