

CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Human IL-17B Kit
Kit Catalog Numbers: K151WO-Series
Kit Lot Number: K0080758
Expiration Date: 28 Feb 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-17B	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-17B	1.12	75% - 125%	< 20%	1040	75% - 125%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IL-17B	Calibrator	1400	pg/mL
		Cytokine Panel 2 (Human) Control 1	414	pg/mL
		Cytokine Panel 2 (Human) Control 2	83.6	pg/mL
		Cytokine Panel 2 (Human) Control 3	30.6	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu IL-17B - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1400	241,000	5.0
350	60,000	3.2
87.5	14,000	8.8
21.9	3,710	5.2
5.47	1,010	2.2
1.37	301	2.7
0.342	113	3.9
0.00	56	5.2



CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-17B Plate	Z0046221	2-8°C	28 Feb 2019
Anti-hu IL-17B Antibody	D0080767	2-8°C	30 Apr 2019
Cytokine Panel 2 (human) Calibrator Blend	A0080116	2-8°C	31 Jan 2022
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), 1000mL	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		21 JUL 2017