

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-27 Kit
Kit Catalog Numbers: K151WX-Series
Kit Lot Number: K0081485
Expiration Date: 31 May 2022

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	IL-27	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-27	38.7	80% - 120%	< 20%	13000	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-27	Calibrator	19300	pg/mL
		TH17 Panel 1 (human) Control 1	31000	pg/mL
		TH17 Panel 1 (human) Control 2	6560	pg/mL
		TH17 Panel 1 (human) Control 3	1380	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IL-27 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
19300	100,000	1.7
4830	28,900	3.2
1210	7,750	4.3
302	2,010	5.6
75.4	604	4.8
18.8	229	4.2
4.71	137	2.8
0.00	111	8.5

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-27 Plate	Z0047324	2-8°C	31 Oct 2022
Anti-hu IL-27 Antibody	D0081429	2-8°C	31 Jul 2023
TH17 Panel 1 (human) Calibrator Blend	A0080219	2-8°C	30 Apr 2024
Diluent 43	M0430036	≤ -10°C	31 May 2022
Diluent 3	M0030096	≤ -10°C	30 Jun 2022
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	Karen Hamilla	Quality		11 JUN 2020