

**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:** K0081449  
**Expiration Date:** 28 Feb 2021

**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	21000	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
21000	47,900	3.4
5250	14,700	4.7
1310	3,970	3.0
328	1,070	2.8
82.0	353	3.7
20.5	159	4.2
5.13	112	6.9
0.00	94	4.3

## CERTIFICATE OF ANALYSIS

Page 2 of 2


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-27 Plate	Z0047203	2-8°C	31 May 2022
Anti-hu IL-27 Antibody	D0081429	2-8°C	31 Jul 2023
TH17 Panel 1 (human) Calibrator Blend	A0080232	2-8°C	31 Oct 2024
Diluent 43	M0430029	≤ -10°C	30 Apr 2021
Diluent 3	M0030085	≤ -10°C	28 Feb 2021
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		13 APR 2020