

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: K0081201
Expiration Date: 30 Jun 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	16100	pg/mL
		TH17 Panel 1 (human) Control 1	18400	pg/mL
		TH17 Panel 1 (human) Control 2	3450	pg/mL
		TH17 Panel 1 (human) Control 3	790	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
16100	64,800	2.7
4030	12,800	6.9
1010	3,160	3.1
252	803	3.3
62.9	258	5.6
15.7	114	4.6
3.93	84	9.3
0.00	73	5.0

CERTIFICATE OF ANALYSIS

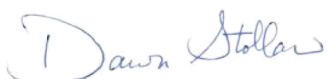
Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-27 Plate	Z0047038	2-8°C	31 Aug 2021
Anti-hu IL-27 Antibody	D0081139	2-8°C	28 Feb 2022
TH17 Panel 1 (human) Calibrator Blend	A0080119	2-8°C	31 Jan 2022
Diluent 43	M0430030	≤ -10°C	30 Jun 2021
Diluent 3	M0030087	≤ -10°C	30 Jun 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), 100 mL	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	Dawn Stollar	Quality		22 APR 2019