

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: K0081525
Expiration Date: 30 Jun 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	21000	pg/mL
		TH17 Panel 1 (human) Control 1	17500	pg/mL
		TH17 Panel 1 (human) Control 2	4200	pg/mL
		TH17 Panel 1 (human) Control 3	826	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	89,400	1.8
5250	25,000	2.6
1310	6,340	1.7
328	1,680	1.3
82.0	482	1.4
20.5	177	2.3
5.13	103	5.0
0.00	79	4.6

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-27 Plate	Z0047368	2-8°C	31 Jan 2023
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	M0430037	≤ -10°C	31 Aug 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), 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	Karen Hamilla	Quality		09 SEP 2020