

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-17A Kit

Kit Catalog Numbers: K152RF-Series

Kit Lot Number: K0081448

Expiration Date: 31 Mar 2022

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		TH17 Panel 1 (mouse) Control 1		TH17 Panel 1 (mouse) Control 2		TH17 Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-17A	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-17A	0.255	75% - 125%	< 20%	360	75% - 125%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-17A	Calibrator	559	pg/mL
		TH17 Panel 1 (mouse) Control 1	508	pg/mL
		TH17 Panel 1 (mouse) Control 2	85.2	pg/mL
		TH17 Panel 1 (mouse) Control 3	14.8	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IL-17A - Spot 1		
Conc. pg/mL	ECL Signal	%CV
559	190,000	2.1
140	49,600	2.6
34.9	12,600	2.4
8.73	3,210	3.1
2.18	839	2.5
0.546	253	3.2
0.136	106	4.1
0.00	55	9.0

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-17A Plate	Z0047280	2-8°C	31 Jul 2022
Anti-ms IL-17A Antibody	D0081190	2-8°C	31 Jul 2022
TH17 Panel 1 (mouse) Calibrator Blend	A0080220	2-8°C	30 Jun 2024
Diluent 41	M0410030	≤ -10°C	31 Mar 2022
Diluent 45	M0450031	≤ -10°C	31 Mar 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 (mouse) Control 1	Not Kit Specific
TH17 Panel 1 (mouse) Control 2	Not Kit Specific
TH17 Panel 1 (mouse) 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		20 MAR 2020