



RFE MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX Plus Human IL-17A Gen. B Kit

Catalog Number(s) / Size(s): K151WM-Series

Lot Number: K0081437

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0081437 has been extended to 31 JUL 2020.

The expiration dates for the following kit components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Human IL-17A Gen. B Plate	Z0046395	31 JUL 2019	31 JUL 2020

Function	Name	Signature	Date
Quality	Karen Hamilla		01 APR 2020

This memo supersedes all other statements regarding date of expiration for this product.

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-17A Gen. B Kit
Kit Catalog Numbers: K151WM-Series
Kit Lot Number: K0081437
Expiration Date: 31 Jul 2019

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-17A Gen. B	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-17A Gen. B	5.86	80% - 120%	< 20%	1950	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-17A Gen. B	Calibrator	2570	pg/mL
		TH17 Panel 1 (human) Control 1	1980	pg/mL
		TH17 Panel 1 (human) Control 2	293	pg/mL
		TH17 Panel 1 (human) Control 3	42.5	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IL-17A Gen. B - Spot 1		
Conc. pg/mL	ECL Signal	%CV
2570	172,000	3.8
643	41,600	2.4
161	9,920	5.6
40.2	2,640	2.7
10.0	717	2.6
2.51	240	2.1
0.627	119	4.3
0.00	80	5.3

CERTIFICATE OF ANALYSIS

Components:


Description	Lot Number	Storage Temperature	Expiration Date
Human IL-17A Gen. B Plate	Z0046395	2-8°C	31 Jul 2019
Anti-hu IL-17A Antibody	D0081135	2-8°C	30 Apr 2022
TH17 Panel 1 (human) Calibrator Blend	A0080232	2-8°C	31 Oct 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), 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		01 APR 2020