



RFE MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX Plus Human IL-17A/F Kit

Catalog Number(s) / Size(s): K151WN-Series

Lot Number: K0080624

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0080624 has been extended to 28 Feb 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/F Plate	Z0046220	28 Feb 2019	28 Feb 2020
Anti-hu IL-17 A/F Antibody	D0080766	30 Apr 2019	28 Feb 2020
Diluent 43	M0430020	30 Jun 2019	28 Feb 2020
Diluent 3	M0030074	30 Jun 2019	28 Feb 2020

Function	Name	Signature	Date
Quality Assurance	Dawn Stollar	<i>Dawn Stollar</i>	14 DEC 2018

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/F Kit
Kit Catalog Numbers: K151WN-Series
Kit Lot Number: K0080624
Expiration Date: 28 Feb 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Cytokine Panel 2 (human) Control 1		Cytokine Panel 2 (human) Control 2		Cytokine Panel 2 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu IL-17A/F	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu IL-17A/F	7.57	80% - 120%	< 20%	3900	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IL-17A/F	Calibrator	5430	pg/mL
		Cytokine Panel 2 (human) Control 1	2390	pg/mL
		Cytokine Panel 2 (human) Control 2	317	pg/mL
		Cytokine Panel 2 (human) Control 3	39.2	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu IL-17A/F - Spot 1		
Conc. pg/mL	ECL Signal	%CV
5430	436,000	1.1
1360	108,000	1.3
339	25,800	2.2
84.8	6,190	1.0
21.2	1,630	1.8
5.30	467	1.9
1.33	178	3.4
0.00	76	5.0

CERTIFICATE OF ANALYSIS

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-17A/F Plate	Z0046220	2-8°C	28 Feb 2019
Anti-hu IL-17A/F antibody	D0080766	2-8°C	30 Apr 2019
Cytokine panel 2 (human) Calibrator Blend	A0080114	2-8°C	31 Dec 2021
Diluent 43	M0430020	≤ -10°C	30 Jun 2019
Diluent 3	M0030074	≤ -10°C	30 Jun 2019
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
Cytokine Panel 2 (human) Control 1	Not Kit Specific
Cytokine Panel 2 (human) Control 2	Not Kit Specific
Cytokine Panel 2 (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 Assurance	<i>Dawn Stollar</i>	21 JUL 2017