



RFE MEMO

Page 1 of 1

Product Description(s): V-Plex® and V-Plex Plus Human IL-17C Kit


Catalog Number(s) / Size(s): K151WP-Series

Lot Number: K0081229

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0081229 has been extended to 31 MAY 2020.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Human IL-17C Plate	Z0046316	31 MAY 2019	31 MAY 2020

Function	Name	Signature	Date
Quality	Karen Hamilla		21 JUN 2019

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-17C Kit
Kit Catalog Numbers: K151WP-Series
Kit Lot Number: K0081229
Expiration Date: 31 May 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	IL-17C	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-17C	3.84	75% - 125%	< 20%	1620	75% - 125%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-17C	Calibrator	2500	pg/mL
		Cytokine Panel 2 (human) Control 1	2120	pg/mL
		Cytokine Panel 2 (human) Control 2	154	pg/mL
		Cytokine Panel 2 (human) Control 3	78.6	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IL-17C - Spot 1		
Conc. pg/mL	ECL Signal	%CV
2500	120,000	4.2
625	34,500	6.4
156	8,200	4.6
39.1	2,100	4.9
9.77	541	3.5
2.44	194	4.1
0.610	102	5.6
0.00	68	5.4

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-17C Plate	Z0046316	2-8°C	31 May 2019
Anti-hu IL-17C Antibody	D0081015	2-8°C	30 Jun 2020
Cytokine Panel 2 (human) Calibrator Blend	A0080116	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), 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	Karen Hamilla	Quality		21 JUN 2019