



**RFE MEMO**

Page 1 of 1

**Product Description(s):** V-PLEX® and V-PLEX Plus Human IL-15 Kit


**Catalog Number(s) / Size(s):** K151RD-Series, K156RD-Series

**Lot Number:** K0081331

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0081331 been extended to 30 SEP 2022.

The expiration date for the following kit component has been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Human IL-15 Plate	Z0047181	28 FEB 2022	30 SEP 2022
Anti-hu IL-15 Antibody	D0081081	30 SEP 2021	30 SEP 2022
Diluent 43	M0430034	28 FEB 2022	30 SEP 2022
Diluent 3	M0030092	31 MAR 2022	30 SEP 2022

Function	Name	Signature	Date
Quality	Karen Hamilla		14 APR 2022

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-15 Kit  
**Kit Catalog Numbers:** K151RD-Series, K156RD-Series  
**Kit Lot Number:** K0081331  
**Expiration Date:** 30 Sep 2021

### Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Cytokine Panel 1 (human) Control 1		Cytokine Panel 1 (human) Control 2		Cytokine Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-15	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-15	0.774	80% - 120%	< 20%	525	80% - 120%	< 20%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-15	Calibrator	976	pg/mL
		Cytokine Panel 1 (human) Control 1	259	pg/mL
		Cytokine Panel 1 (human) Control 2	40.7	pg/mL
		Cytokine Panel 1 (human) Control 3	7.82	pg/mL

Controls tested may not be the same lot as shipped.

### Standard Curve Data:

IL-15 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
976	222,000	3.7
244	39,200	3.2
61.0	8,850	3.2
15.3	2,310	2.4
3.81	617	3.0
0.953	204	3.1
0.238	95	3.7
0.00	61	5.9

## CERTIFICATE OF ANALYSIS

Page 2 of 2


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-15 Plate	Z0047181	2-8°C	28 Feb 2022
Anti-hu IL-15 Antibody	D0081081	2-8°C	30 Sep 2021
Cytokine Panel 1 (human) Calibrator Blend	A0080227	2-8°C	31 Oct 2024
Diluent 43	M0430034	≤ -10°C	28 Feb 2022
Diluent 3	M0030092	≤ -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
Cytokine Panel 1 (human) Control 1	Not Kit Specific
Cytokine Panel 1 (human) Control 2	Not Kit Specific
Cytokine 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
<b>Review/Approval</b>	Karen Hamilla	Quality		29 JAN 2020