



RFE MEMO

Page 1 of 1

Product Description(s): V-Plex® and V-Plex Plus Mouse IL-33 Kit

Catalog Number(s) / Size(s): K152XB-Series

Lot Number: K0080742

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0080742 has been extended to 31 Aug 2020.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Mouse IL-33 Plate	Z0046398	31 Aug 2019	31 Aug 2020
Anti-ms IL-33 Antibody	D0080938	31 Aug 2019	31 Aug 2020
Diluent 41	M0410016	31 Aug 2019	31 Aug 2020
Diluent 45	M0450016	30 Sep 2019	31 Aug 2020

Function	Name	Signature	Date
Quality	Karen Hamilla		14 OCT 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 Mouse IL-33 Kit

Kit Catalog Numbers: K152XB-Series

Kit Lot Number: K0080742

Expiration Date: 31 Aug 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Cytokine Panel 1 (mouse) Control 1		Cytokine Panel 1 (mouse) Control 2		Cytokine Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	ms IL-33	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	ms IL-33	1.85	75% - 125%	< 20%	1950	75% - 125%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms IL-33	Calibrator	3280	pg/mL
		Cytokine Panel 1 (mouse) Control 1	2450	pg/mL
		Cytokine Panel 1 (mouse) Control 2	290	pg/mL
		Cytokine Panel 1 (mouse) Control 3	36.8	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

ms IL-33 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
3280	518,000	3.0
820	153,000	2.4
205	40,800	3.3
51.3	10,100	2.0
12.8	2,740	2.2
3.20	770	3.1
0.801	294	3.7
0.00	126	3.3

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-33 Plate	Z0046398	2-8°C	31 Aug 2019
Anti-ms IL-33 Antibody	D0080938	2-8°C	31 Aug 2019
Cytokine Panel 1 (mouse) Kit Calibrator Blend	A0080125	2-8°C	31 Mar 2022
Diluent 41	M0410016	≤ -10°C	31 Aug 2019
Diluent 45	M0450016	≤ -10°C	30 Sep 2019
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
MSD Wash Buffer (20x), 100 mL	Not Kit Specific
Plate Seals	Not Kit Specific
Cytokine Panel 1 (mouse) Control 1	Not Kit Specific
Cytokine Panel 1 (mouse) Control 2	Not Kit Specific
Cytokine 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	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	28 JUL 2017