



RFE MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX Plus Mouse IFN- γ Kit

Catalog Number(s) / Size(s): K152QO-Series

Lot Number: K0080974

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0080974 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
Anti-ms IFN- γ Antibody	D0080851	31 May 2019	31 May 2020

Function	Name	Signature	Date
Quality	Karen Hamilla		25 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 Mouse IFN- γ Kit
Kit Catalog Numbers: K152QO-Series
Kit Lot Number: K0080974
Expiration Date: 31 May 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (mouse) Control 1		Proinflammatory Panel 1 (mouse) Control 2		Proinflammatory Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Mouse IFN- γ	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Mouse IFN- γ	0.390	75% - 125%	< 25%	570	75% - 125%	< 25%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Mouse IFN- γ	Calibrator	938	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	282	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	21.5	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	1.76	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

Mouse IFN- γ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
938	441,000	3.1
235	102,000	3.7
58.6	24,500	2.1
14.7	5,950	4.3
3.66	1,670	2.4
0.916	496	2.9
0.229	197	3.4
0.00	96	5.9

CERTIFICATE OF ANALYSIS

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IFN- γ Plate	Z0046781	2-8°C	31 Aug 2020
Anti-ms IFN- γ Antibody	D0080851	2-8°C	31 May 2019
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080138	2-8°C	31 May 2022
Diluent 41	M0410022	$\leq -10^{\circ}\text{C}$	31 Jul 2020
Diluent 45	M0450020	$\leq -10^{\circ}\text{C}$	31 May 2020
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
Proinflammatory Panel 1 (mouse) Control 1	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 2	Not Kit Specific
Proinflammatory 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>	11 MAY 2018