



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: K0081049

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0081049 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
Anti-ms IFN- γ Antibody	D0080960	31 AUG 2019	31 AUG 2020
Diluent 41	M0410022	31 JUL 2020	31 AUG 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

Product Description: V-PLEX® and V-PLEX Plus Mouse IFN- γ Kit
Kit Catalog Numbers: K152QO-Series
Kit Lot Number: K0081049
Expiration Date: 31 Aug 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	ms 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	ms IFN- γ	0.390	75% - 125%	< 25%	570	75% - 125%	< 25%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms IFN- γ	Calibrator	644	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	467	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	32.2	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	2.15	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

ms IFN- γ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
644	303,000	13.0
161	74,100	6.1
40.3	17,600	6.9
10.1	4,240	6.5
2.52	1,120	4.8
0.629	329	5.9
0.157	135	5.5
0.00	71	5.4

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IFN- γ Plate	Z0046869	2-8°C	31 Dec 2020
Anti-ms IFN- γ Antibody	D0080960	2-8°C	31 Aug 2019
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080152	2-8°C	30 Nov 2022
Diluent 41	M0410022	\leq -10°C	31 Jul 2020
Diluent 45	M0450021	\leq -10°C	30 Sep 2020
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
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		03 OCT 2018