



**RFE MEMO**

Page 1 of 1

**Product Description(s):** V-PLEX® and V-PLEX Plus Mouse IL-12p70 Kit

**Catalog Number(s) / Size(s):** K152QV - Series

**Lot Number:** K0080849

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0080849 has been extended to 30 APR 2019.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Mouse IL-12p70 Plate	Z0045953	30 APR 2018	30 APR 2019

Function	Name	Signature	Date
Quality Assurance	Dawn Stollar	<i>Dawn Stollar</i>	12 MAR 2018

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-12p70 Kit  
**Kit Catalog Numbers:** K152QV-Series  
**Kit Lot Number:** K0080849  
**Expiration Date:** 30 Apr 2018

**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 IL-12p70	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	ms IL-12p70	179	75% - 125%	< 25%	20600	75% - 125%	< 25%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	ms IL-12p70	Calibrator	31500	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	13400	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	3840	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	1120	pg/mL

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

ms IL-12p70 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
31500	508,000	2.0
7880	95,600	4.0
1970	20,700	3.0
492	4,970	2.9
123	1,400	3.7
30.8	428	2.3
7.69	205	3.3
0.00	94	5.4

**CERTIFICATE OF ANALYSIS**

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-12p70 Plate	Z0045953	2-8°C	30 Apr 2018
Anti-ms IL-12p70 Antibody	D0080879	2-8°C	30 Jun 2019
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080110	2-8°C	31 Dec 2021
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	
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
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	10 AUG 2017