



**RFE MEMO**

Page 1 of 1

**Product Description(s):** V-PLEX® and V-PLEX Plus Mouse KC/GRO Kit

**Catalog Number(s) / Size(s):** K152QT-Series

**Lot Number:** K0080705

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0080705 has been extended to 31 Mar 2019.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Diluent 45	M0450012	28 FEB 2019	31 MAR 2019
Anti-ms KC/GRO Antibody	D0080639	31 MAR 2018	31 MAR 2019

Function	Name	Signature	Date
Quality Assurance	Dawn Stollar	<i>Dawn Stollar</i>	15 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 KC/GRO Kit  
**Kit Catalog Numbers:** K152QT-Series  
**Kit Lot Number:** K0080705  
**Expiration Date:** 31 Mar 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 KC/GRO	80% - 120%	< 20%	80% - 130%	< 25%	80% - 130%	< 25%	80% - 130%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	ms KC/GRO	3.29	75% - 125%	< 25%	1230	75% - 125%	< 25%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	ms KC/GRO	Calibrator	2110	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	902	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	111	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	16.2	pg/mL

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

ms KC/GRO - Spot 1		
Conc. pg/mL	ECL Signal	%CV
2110	660,000	1.8
528	141,000	1.6
132	31,400	2.1
33.0	7,540	1.9
8.24	1,940	1.8
2.06	559	2.4
0.515	204	2.0
0.00	85	3.7

**CERTIFICATE OF ANALYSIS**


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Mouse KC/GRO Plate	Z0046297	2-8°C	30 Jun 2019
Anti-ms KC/GRO Antibody	D0080639	2-8°C	31 Mar 2018
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080071	2-8°C	30 Nov 2020
Diluent 41	M0410015	≤ -10°C	30 Jun 2019
Diluent 45	M0450012	≤ -10°C	28 Feb 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)	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		15 FEB 2017