



RFE MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX Plus Human MCP-1 Kit

Catalog Number(s) / Size(s): K151NN-Series

Lot Number: K0080876

Based on additional testing, the original date of expiration for this kit has been extended by 6 months. The expiration date of K0080876 has been extended to 31 Dec 2019.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Anti-hu MCP-1 Antibody	D0080865	30 JUN 2019	31 DEC 2019

Function	Name	Signature	Date
Quality Assurance	Dawn Stollar	<i>Dawn Stollar</i>	12 FEB 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 Human MCP-1 Kit

Kit Catalog Numbers: K151NN-Series

Kit Lot Number: K0080876

Expiration Date: 30 Jun 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Chemokine Panel 1(Human) Control 1		Chemokine Panel 1(Human) Control 2		Chemokine Panel 1(Human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu MCP-1	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu MCP-1	1.09	80% - 120%	< 20%	375	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu MCP-1	Calibrator	686	pg/mL
		Chemokine Panel 1(Human) Control 1	420	pg/mL
		Chemokine Panel 1(Human) Control 2	66.6	pg/mL
		Chemokine Panel 1(Human) Control 3	11.1	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu MCP-1 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
686	212,000	3.6
172	53,300	3.3
42.9	13,500	2.0
10.7	3,490	2.8
2.68	943	3.4
0.670	277	2.5
0.167	110	2.1
0.00	56	5.6

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MCP-1 Plate	Z0046607	2-8°C	31 Jan 2020
Anti-hu MCP-1 Antibody	D0080865	2-8°C	30 Jun 2019
Chemokine Panel 1 (Human) Calibrator Blend	A0080131	2-8°C	31 Jul 2022
Diluent 43	M0430023	≤ -10°C	31 Dec 2019
Diluent 3	M0030078	≤ -10°C	31 Jan 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
Chemokine Panel 1(Human) Control 1	Not Kit Specific
Chemokine Panel 1(Human) Control 2	Not Kit Specific
Chemokine Panel 1(Human) 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		10 OCT 2017