



**RFE MEMO**

Page 1 of 1

**Product Description(s):** V-PLEX® and V-PLEX Plus NHP MIP-1β Kit

**Catalog Number(s) / Size(s):** K156NR-Series

**Lot Number:** K0081062

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0081062 has been extended to 30 SEP 2019.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Human MIP-1β Plate	Z0046116	30 SEP 2018	30 SEP 2019
Anti-hu MIP-1β Antibody	D0080861	30 JUN 2019	30 SEP 2019

Function	Name	Signature	Date
Quality Assurance	Dawn Stollar	<i>Dawn Stollar</i>	10 OCT 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 NHP MIP-1 $\beta$  Kit  
**Kit Catalog Numbers:** K156NR-series  
**Kit Lot Number:** K0081062  
**Expiration Date:** 30 SEP 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	NHP MIP-1 $\beta$	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	NHP MIP-1 $\beta$	1.02	80% - 120%	< 20%	750	80% - 120%	< 20%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	NHP MIP-1 $\beta$	Calibrator	1190	pg/mL
		Chemokine Panel 1 (human) Control 1	1000	pg/mL
		Chemokine Panel 1 (human) Control 2	176	pg/mL
		Chemokine Panel 1 (human) Control 3	25.8	pg/mL

Controls tested may not be the same lot as shipped.

The calibrator and controls for this assay contain a human recombinant cytokine calibrator sequence that is highly homologous to the non-human primate cytokine. This assay has been verified to cross-react with homologs of rhesus and cynomolgus monkeys. See Product Insert for experimental data.

### Standard Curve Data:

NHP MIP-1 $\beta$ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1190	642,000	2.4
298	152,000	2.2
74.4	25,500	1.5
18.6	4,220	2.6
4.65	901	1.9
1.16	267	2.7
0.291	123	3.0
0.00	78	5.6

## CERTIFICATE OF ANALYSIS

Page 2 of 2

### Components:


Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 $\beta$ Plate	Z0046116	2-8°C	30 Sep 2019
Anti-hu MIP-1 $\beta$ Antibody	D0080861	2-8°C	30 Jun 2019
Chemokine Panel 1 (human) Calibrator Blend	A0080079	2-8°C	31 May 2021
Diluent 43	M0430021	$\leq$ -10°C	30 Sep 2019
Diluent 3	M0030077	$\leq$ -10°C	30 Nov 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
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
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		10 OCT 2018