



**RFE MEMO**

Page 1 of 1

**Product Description(s):** V-PLEX® and V-PLEX Plus Human IP-10 Kit

**Catalog Number(s) / Size(s):** K151NV-Series

**Lot Number:** K0081018

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0081018 has been extended to 31 Dec 2020.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Human IP-10 Plate	Z0046839	30 NOV 2020	31 DEC 2020
Anti-hu IP-10 Antibody	D0080994	31 DEC 2019	31 DEC 2020
Diluent 43	M0430026	31 AUG 2020	31 DEC 2020
Diluent 3	M0030082	31 OCT 2020	31 DEC 2020

Function	Name	Signature	Date
Quality	Karen Hamilla		18 OCT 2019

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 IP-10 Kit  
**Kit Catalog Numbers:** K151NV-Series  
**Kit Lot Number:** K0081018  
**Expiration Date:** 31 Dec 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 IP-10	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu IP-10	1.37	80% - 120%	< 20%	500	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	hu IP-10	Calibrator	2090	pg/mL
		Chemokine Panel 1 (human) Control 1	601	pg/mL
		Chemokine Panel 1 (human) Control 2	92.1	pg/mL
		Chemokine Panel 1 (human) Control 3	15.5	pg/mL

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

hu IP-10 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
2090	418,000	2.2
523	126,000	1.9
131	32,400	3.6
32.7	8,360	2.8
8.16	2,130	3.4
2.04	600	2.4
0.510	189	2.1
0.00	66	5.4

## CERTIFICATE OF ANALYSIS

Page 2 of 2


### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IP-10 Plate	Z0046839	2-8°C	30 Nov 2020
Anti-hu IP-10 Antibody	D0080994	2-8°C	31 Dec 2019
Chemokine Panel 1 (human) Calibrator Blend	A0080171	2-8°C	30 Jun 2023
Diluent 43	M0430026	≤ -10°C	31 Aug 2020
Diluent 3	M0030082	≤ -10°C	31 Oct 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), 100 mL	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 SEP 2018