



RFE MEMO

Page 1 of 1

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

Catalog Number(s) / Size(s): K151RH-Series, K156RH-Series

Lot Number: K0080879

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0080879 has been extended to 31 JUL 2020.

The expiration date for the following kit component has been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Human VEGF Plate	Z0056036	31 JAN 2020	31 JUL 2020
Anti-hu VEGF Antibody	D0080894	31 JUL 2019	31 JUL 2020
Diluent 43	M0430023	31 DEC 2019	31 JUL 2020
Diluent 3	M0030078	31 JAN 2020	31 JUL 2020

Function	Name	Signature	Date
Quality	Karen Hamilla		21 JAN 2020

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 VEGF Kit
Kit Catalog Numbers: K156RH-Series
Kit Lot Number: K0080879
Expiration Date: 31 Jul 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Cytokine Panel 1 (human) Control 1		Cytokine Panel 1 (human) Control 2		Cytokine Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	NHP VEGF	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	NHP VEGF	7.70	80% - 120%	< 20%	562	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	NHP VEGF	Calibrator	1340	pg/mL
		Cytokine Panel 1 (human) Control 1	373	pg/mL
		Cytokine Panel 1 (human) Control 2	127	pg/mL
		Cytokine Panel 1 (human) Control 3	56.1	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 VEGF - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1340	286,000	1.3
335	56,300	1.8
83.8	11,800	1.9
20.9	2,720	1.0
5.23	729	2.7
1.31	220	3.1
0.327	95	3.4
0.00	53	8.1

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human VEGF Plate	Z0056036	2-8°C	31 Jan 2020
Anti-hu VEGF Antibody	D0080894	2-8°C	31 Jul 2019
Cytokine Panel 1 (human) Calibrator Blend	A0080108	2-8°C	30 Nov 2021
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), 100 mL	Not Kit Specific
Cytokine Panel 1 (human) Control 1	Not Kit Specific
Cytokine Panel 1 (human) Control 2	Not Kit Specific
Cytokine 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	<i>Dawn Stollar</i>	01 DEC 2017