



RFE MEMO

Page 1 of 1

Product Description(s): V-PLEX© and V-PLEX Plus NHP Eotaxin-3 Kit


Catalog Number(s) / Size(s): K156NU-Series

Lot Number: K0080748

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0080748 has been extended to 30 JUN 2020.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Human Eotaxin-3 Plate	Z0046423	31 AUG 2019	30 JUN 2020
Diluent 43	M0420020	30 JUN 2019	30 JUN 2020
Diluent 3	M0030074	30 JUN 2019	30 JUN 2020
Anti-hu Eotaxin-3 Antibody	D0080862	30 JUN 2019	30 JUN 2020

Function	Name	Signature	Date
Quality Assurance	Dawn Stollar		01 APR 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 NHP Eotaxin-3 Kit
Kit Catalog Numbers: K156NU-Series
Kit Lot Number: K0080748
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	NHP Eotaxin-3	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	NHP Eotaxin-3	10.2	80% - 120%	< 20%	3750	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	NHP Eotaxin-3	Calibrator	7120	pg/mL
		Chemokine Panel 1 (human) Control 1	4670	pg/mL
		Chemokine Panel 1 (human) Control 2	794	pg/mL
		Chemokine Panel 1 (human) Control 3	127	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 Eotaxin-3 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
7120	85,000	3.6
1780	27,000	1.4
445	7,750	2.2
111	1,990	2.3
27.8	564	2.8
6.95	188	3.4
1.74	96	5.2
0.00	62	7.9

CERTIFICATE OF ANALYSIS

Page 2 of 2

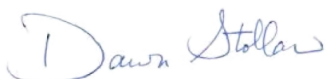
Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human Eotaxin-3 Plate	Z0046423	2-8°C	31 Aug 2019
Anti-hu Eotaxin-3 Antibody	D0080862	2-8°C	30 Jun 2019
Chemokine Panel 1 (human) Calibrator Blend	A0080122	2-8°C	28 Feb 2022
Diluent 43	M0430020	≤ -10°C	30 Jun 2019
Diluent 3	M0030074	≤ -10°C	30 Jun 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
Review/Approval	Dawn Stollar	Quality Assurance		01 APR 2019