



RFE MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-Plex Plus Rat IL-10 Kit

Catalog Number(s) / Size(s): K153QU-series

Lot Number: K0080737

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0080737 has been extended to 30 APR 2019.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Anti-rat IL-10 Antibody	D0080727	31 Oct 2018	30 APR 2019
Blocker H	R0100011	30 Apr 2018	30 APR 2019
Diluent 42	M0420009	31 Aug 2018	30 APR 2019

Function	Name	Signature	Date
Quality Assurance	Dawn Stollar	<i>Dawn Stollar</i>	18 SEP 2018

This memo supersedes all other statements regarding date of expiration for this product.



CERTIFICATE OF ANALYSIS

Page 1 of 3

Product Description: V-PLEX® and V-PLEX Plus Rat IL-10 Kit
Kit Catalog Numbers: K153QU-series
Kit Lot Number: K0080737
Expiration Date: 30 Apr 2018

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (rat) Control 1		Proinflammatory Panel 1 (rat) Control 2		Proinflammatory Panel 1 (rat) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Rat IL-10	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Rat IL-10	163	75% - 125%	< 25%	15700	75% - 125%	< 25%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Rat IL-10	Calibrator	26100	pg/mL
		Proinflammatory Panel 1 (rat) Control 1	19800	pg/mL
		Proinflammatory Panel 1 (rat) Control 2	4360	pg/mL
		Proinflammatory Panel 1 (rat) Control 3	1640	pg/mL

Controls tested may not be the same lot as shipped.



CERTIFICATE OF ANALYSIS

Page 2 of 3

Standard Curve Data:

Rat IL-10 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
26100	908,000	7.5
6530	237,000	5.2
1630	56,600	3.9
408	14,100	4.0
102	3,580	3.5
25.5	968	7.2
6.37	324	3.6
0.00	116	4.9

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Rat IL-10 Plate	Z0046406	2-8°C	31 Aug 2019
Anti-rat IL-10 Antibody	D0080727	2-8°C	31 Oct 2018
Proinflammatory Panel 1 (rat) Calibrator Blend	A0080083	2-8°C	31 May 2021
Blocker H	R0100011	Room Temperature	30 Apr 2018
Diluent 42	M0420009	≤ -10°C	31 Aug 2018
Diluent 40	M0400010	≤ -10°C	31 May 2019
MSD Read Buffer (4x)	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
Proinflammatory Panel 1 (rat) Control 1	Not Kit Specific
Proinflammatory Panel 1 (rat) Control 2	Not Kit Specific
Proinflammatory Panel 1 (rat) Control 3	Not Kit Specific

Statement:

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.



CERTIFICATE OF ANALYSIS

Page 3 of 3

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	04 APR 2017