



RFE MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX Plus Human IL-10 Kit
Catalog Number(s) / Size(s): K151QU-Series
Lot Number: K0080354

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0080354 has been extended to 31 Jan 2018.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Human IL-10 Plate	Z0045754	30 Nov 2017	31 Jan 2018
Anti-hu IL-10 Antibody	D0080531	30 Sep 2017	31 Jan 2018
Proinflammatory Panel 1 (human) Calibrator Blend	A0080058	31 May 2020	31 May 2020
Diluent 2	M0020085	31 Jan 2017	31 Jan 2018
Diluent 3	M0030066	30 Sep 2017	31 Jan 2018

Function	Name	Signature	Date
Quality Assurance	Dawn Stollar	<i>Dawn Stollar</i>	08 FEB 2017

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



CERTIFICATE OF ANALYSIS

COA-5717

Page 1 of 3

Product Description: Human IL-10 Kit V-PLEX® and V-PLEX Plus
Kit Catalog Numbers: K151QUD-1, K151QUD-2, K151QUD-4, K151QUG-1, K151QUG-2, K151QUG-4
Kit Lot Number: K0080354
Expiration Date: 31 Jan 2017

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (human) Control 1		Proinflammatory Panel 1 (human) Control 2		Proinflammatory Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu IL-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 IL-10	0.680	80% - 120%	< 20%	233	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IL-10	Calibrator	317	pg/mL
		Proinflammatory Panel 1 (human) Control 1	95.8	pg/mL
		Proinflammatory Panel 1 (human) Control 2	24.7	pg/mL
		Proinflammatory Panel 1 (human) Control 3	3.69	pg/mL

Controls tested may not be the same lot as shipped.



CERTIFICATE OF ANALYSIS

COA-5717

Page 2 of 3

Standard Curve Data:

hu IL-10 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
317	126,192	1.9
79.3	33,463	1.7
19.8	8,516	1.9
4.95	2,157	2.2
1.24	577	2.3
0.310	192	1.9
0.0774	98	3.2
0	65	5.9

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-10 Plate	Z0045754	2-8°C	30 Nov 2017
Anti-hu IL-10 Antibody	D0080531	2-8°C	30 Sep 2017
Proinflammatory Panel 1 (human) Calibrator Blend	A0080058	2-8°C	31 May 2020
Diluent 2	M0020085	≤ -10°C	31 Jan 2017
Diluent 3	M0030066	≤ -10°C	30 Sep 2017
MSD Read Buffer T (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 (human) Control 1	Not Kit Specific
Proinflammatory Panel 1 (human) Control 2	Not Kit Specific
Proinflammatory Panel 1 (human) Control 3	Not Kit Specific

Statement:


The above kit was tested using MSD procedures and it conforms to MSD's specifications. The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.



CERTIFICATE OF ANALYSIS

COA-5717

Page 3 of 3

	Name	Function	Signature	Date
Review/Approval	Anastasia Belilovets	Quality Assurance		15 Jul 2015