



RFE MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX Plus Human IL-6 Kit

Catalog Number(s) / Size(s): K151QX-Series; K156QX-Series

Lot Number: K0081460

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0081460 has been extended to 31 JUL 2023.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Human IL-6 Plate	Z0047300	31 AUG 2022	31 JUL 2023
Anti-hu IL-6 Antibody	D0081311	30 NOV 2022	31 JUL 2023
Diluent 2	M0020112	31 JUL 2022	31 JUL 2023
Diluent 3	M0030097	31 AUG 2022	31 JUL 2023

Function	Name	Signature	Date
Quality	Karen Hamilla		21 SEP 2022

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 IL-6 Kit
Kit Catalog Numbers: K151QX-Series; K156QX-Series
Kit Lot Number: K0081460
Expiration Date: 31 Jul 2022

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	IL-6	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-6	0.633	80% - 120%	< 20%	488	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-6	Calibrator	755	pg/mL
		Proinflammatory Panel 1 (human) Control 1	289	pg/mL
		Proinflammatory Panel 1 (human) Control 2	49.1	pg/mL
		Proinflammatory Panel 1 (human) Control 3	7.89	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IL-6 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
755	369,000	1.2
189	80,700	1.8
47.2	17,600	3.3
11.8	4,250	2.7
2.95	1,110	2.4
0.737	311	2.3
0.184	123	4.7
0.00	71	16.0

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-6 Plate	Z0047300	2-8°C	31 Aug 2022
Anti-hu IL-6 Antibody	D0081311	2-8°C	30 Nov 2022
Proinflammatory Panel 1 (human) Calibrator Blend	A0080216	2-8°C	30 Jun 2024
Diluent 2	M0020112	≤ -10°C	31 Jul 2022
Diluent 3	M0030097	≤ -10°C	31 Aug 2022
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
MSD Wash Buffer (20X), 100mL	Not Kit Specific
Plate Seals	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 product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		20 APR 2020