



RFE MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX Plus Human IL-5 Kit

Catalog Number(s) / Size(s): K151QS-Series

Lot Number: K0081028

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0081028 has been extended to 31 JUL 2020.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Anti-hu IL-5 Antibody	D0080887	31 JUL 2019	31 JUL 2020

Function	Name	Signature	Date
Quality Assurance	Dawn Stollar		29 MAR 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 Human IL-5 Kit

Kit Catalog Numbers: K151QS-Series

Kit Lot Number: K0081028

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	hu IL-5	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-5	4.41	80% - 120%	< 20%	562	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IL-5	Calibrator	844	pg/mL
		Cytokine Panel 1 (human) Control 1	310	pg/mL
		Cytokine Panel 1 (human) Control 2	110	pg/mL
		Cytokine Panel 1 (human) Control 3	34	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu IL-5 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
844	269,000	1.6
211	72,900	2.5
52.8	18,900	2.4
13.2	4,790	2.7
3.30	1,280	1.3
0.824	380	2.1
0.206	147	3.0
0.00	66	6.5

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-5 Plate	Z0046849	2-8°C	30 Nov 2020
Anti-hu IL-5 Antibody	D0080887	2-8°C	31 Jul 2019
Cytokine Panel 1 (human) Calibrator Blend	A0080121	2-8°C	28 Feb 2022
Diluent 43	M0430026	≤ -10°C	31 Aug 2020
Diluent 3	M0030082	≤ -10°C	31 Oct 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), 100mL	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		20 JUL 2018