

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-5 Kit
Kit Catalog Numbers: K151QS-Series; K156QS-Series
Kit Lot Number: K0081283
Expiration Date: 31 Dec 2021

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

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-5	Calibrator	868	pg/mL
		Cytokine Panel 1 (human) Control 1	326	pg/mL
		Cytokine Panel 1 (human) Control 2	103	pg/mL
		Cytokine Panel 1 (human) Control 3	32.2	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IL-5 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
868	149,000	1.8
217	36,500	2.7
54.3	9,180	2.6
13.6	2,360	1.8
3.39	613	4.2
0.848	195	2.8
0.212	88	4.7
0.00	56	6.4

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-5 Plate	Z0047134	2-8°C	31 Dec 2021
Anti-hu IL-5 Antibody	D0081184	2-8°C	30 Apr 2022
Cytokine Panel 1 (human) Calibrator Blend	A0080188	2-8°C	30 Nov 2023
Diluent 43	M0430033	≤ -10°C	31 Jan 2022
Diluent 3	M0030090	≤ -10°C	31 Dec 2021
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	Karen Hamilla	Quality		20 SEP 2019