

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: K0081472

Expiration Date: 31 Aug 2022

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	895	pg/mL
		Cytokine Panel 1 (human) Control 1	334	pg/mL
		Cytokine Panel 1 (human) Control 2	107	pg/mL
		Cytokine Panel 1 (human) Control 3	33.1	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
895	278,000	2.5
224	71,500	3.3
55.9	18,100	2.5
14.0	4,570	2.3
3.50	1,170	3.1
0.874	346	2.4
0.219	135	4.5
0.00	65	6.8

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-5 Plate	Z0047314	2-8°C	30 Sep 2022
Anti-hu IL-5 Antibody	D0081416	2-8°C	30 Apr 2023
Cytokine Panel 1 (human) Calibrator Blend	A0080227	2-8°C	31 Oct 2024
Diluent 43	M0430037	≤ -10°C	31 Aug 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	
Plate Seals	Not Kit Specific
Wash Buffer (20X), 100 mL	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		26 MAY 2020