

## 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:** K0081382  
**Expiration Date:** 31 Dec 2020

### 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	20% - 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	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
895	279,000	3.1
224	70,600	2.2
55.9	17,500	1.9
14.0	4,330	1.9
3.50	1,120	1.4
0.874	319	2.5
0.219	119	2.5
0.00	57	8.0

## CERTIFICATE OF ANALYSIS

Page 2 of 2


### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-5 Plate	Z0047232	2-8°C	31 May 2022
Anti-hu IL-5 Antibody	D0081361	2-8°C	28 Feb 2023
Cytokine Panel 1 (human) Calibrator Blend	A0080227	2-8°C	31 Oct 2024
Diluent 43	M0430027	≤ -10°C	31 Dec 2020
Diluent 3	M0030084	≤ -10°C	31 Jan 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
<b>Review/Approval</b>	Karen Hamilla	Quality		29 JAN 2020