



# CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-5 Kit  
**Kit Lot Number:** K0082327  
**Expiration Date:** 31 Jul 2025  
**Kit Catalog Numbers:** K151QS-Series; K156QS-Series

### Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-5 Plate	Z0047888	2-8°C	31 Dec 2025
Anti-hu IL-5 Antibody	D0081770	2-8°C	31 Jul 2025
Cytokine Panel 1(human) Calibrator Blend	A0080331	2-8°C	31 Jul 2027
Diluent 3	M0030120	≤ -10°C	31 Jul 2025
Diluent 43	M0430053	≤ -10°C	31 Jan 2026

Additional components may be provided to execute the assay. Additional components are not lot-specific.

### Precision and Accuracy 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 Calibrator:

Spot	Name	Conc.	Units
1	IL-5	888	pg/mL

Dilution to Top of Curve = 1X

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Michelle Likiam	Quality		28 Nov 2023