

**CERTIFICATE OF ANALYSIS**

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-10Kit  
**Lot Number:** K0082525  
**Expiration Date:** 28 FEB 2025  
**Catalog Number:** K151QU-Series; K156QU-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-10 Plate	Z0047789	2-8°C	30 APR 2025
Proinflammatory Panel 1(human) Calibrator Blend	A0080339	2-8°C	31 JUL 2027
Anti-hu IL-10 Antibody	D0081975	2-8°C	30 JUN 2026
Diluent 2	M0020123	≤ -10°C	28 FEB 2025
Diluent 3	M0030118	≤ -10°C	31 MAR 2025

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

**Precision, Accuracy, and Sensitivity Specifications:**

Spot	Name	Calibrator		Proinflammatory Panel 1 (human) Control 1		Proinflammatory Panel 1 (human) Control 2		Proinflammatory Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-10	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-10	0.298	80% - 120%	< 20%	233	80% - 120%	< 20%

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	IL-10	386	pg/mL

Dilution to top of Curve: 1X

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		24 SEP 2024