

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IL-10 Kit

Kit Lot Number: K0082420

Expiration Date: 31 Dec 2024

Kit Catalog Numbers: K151QU-Series, K156QU-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-10 Plate	Z0047726	2-8°C	31 Dec 2024
Anti-hu IL-10 Antibody	D0081876	2-8°C	31 Dec 2025
Proinflammatory Panel 1 (human) Calibrator Blend	A0080312	2-8°C	30 Nov 2026
Diluent 2	M0020126	≤ -10°C	31 Aug 2025
Diluent 3	M0030121	≤ -10°C	31 Oct 2025

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

Precision and Accuracy Specifications:

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


		LLOQ			ULOQ		
Spot	Name	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	377	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	Karen Hamilla	Quality		26 Mar 2024