

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Human IL-10 Kit  
**Kit Lot Number:** K0081629  
**Expiration Date:** 31 Oct 2022  
**Kit Catalog Numbers:** K151QUD-Series; K151QUG-Series

**Precision and Accuracy 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%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 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	380	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-10 Plate	Z0047452	2-8°C	30 Jun 2023
Anti-hu IL-10 Antibody	D0081437	2-8°C	30 Jun 2023
Proinflammatory Panel 1 (human) Calibrator Blend	A0080269	2-8°C	30 Sep 2025
Diluent 2	M0020115	≤ -10°C	31 May 2023
Diluent 3	M0030100	≤ -10°C	31 Oct 2022

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

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		07 JUN 2021