

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IL-10 Kit
Kit Lot Number: K0081999
Expiration Date: 31 Oct 2023
Kit Catalog Numbers: K151QU-Series; K156QU-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	347	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-10 Plate	Z0047682	2-8°C	30 Sep 2024
Anti-hu IL-10 Antibody	D0081553	2-8°C	29 Feb 2024
Proinflammatory Panel 1(human) Calibrator Blend	A0080321	2-8°C	28 Feb 2027
Diluent 2	M0020117	≤ -10°C	31 Oct 2023
Diluent 3	M0030108	≤ -10°C	30 Nov 2023

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		03 JUN 2022