



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Human IL-10 Kit**Lot Number:** K0082853**Expiration Date:** 31 AUG 2027**Catalog Number:** K151QU-Series, K156QU-Series**Lot-Specific Components:**

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IL-10 Antibody	D0082217	2-8°C	31 AUG 2027
Human IL-10, 96-Well Small Spot SECTOR Plate	Z0048210	2-8°C	30 JUN 2028
Proinflammatory Panel 1 (human) Calibrator Blend	A0080414	2-8°C	31 JAN 2030
Diluent 2	M0020134	≤ -10°C	30 SEP 2027
Diluent 3	M0030140	≤ -10°C	30 SEP 2027

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

Precision, Accuracy, and Sensitivity 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	373	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
Review/Approval	Karen Hamilla	Quality		23 JAN 2026