



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Human IL-6 Kit, QuickPlex
Lot Number: K0082644
Expiration Date: 31 MAR 2027
Catalog Number: K151QX-Series, K156QX-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IL-6 Antibody	D0082178	2-8°C	30 JUN 2027
Human IL-6, 96-Well Small Spot QuickPlex Ultra Plate	Z0048077	2-8°C	31 JUL 2027
Proinflammatory Panel 1 (human) Calibrator Blend	A0080365	2-8°C	30 APR 2028
Diluent 2	M0020132	≤ -10°C	31 MAR 2027
Diluent 3	M0030136	≤ -10°C	31 MAY 2027

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-6	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-6	0.633	80% - 120%	< 20%	488	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IL-6	725	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		18 MAR 2025