



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Human IL-1 α Kit**Lot Number:** K0082808**Expiration Date:** 31 MAR 2026**Catalog Number:** K151RB-Series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IL-1 α Antibody	D0081948	2-8°C	31 MAR 2026
Cytokine Panel 1 (human) Calibrator Blend	A0080331	2-8°C	31 JUL 2027
Human IL-1 α 96-Well Small Spot SECTOR Plate	Z0047992	2-8°C	30 NOV 2026
Diluent 3	M0030128	$\leq -10^{\circ}\text{C}$	31 OCT 2026
Diluent 43	M0430055	$\leq -10^{\circ}\text{C}$	30 APR 2026

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

Precision, Accuracy, and Sensitivity Specifications:

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

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-1 α	2.85	80% - 120%	< 20%	278	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IL-1 α	422	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		27 OCT 2025