



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Human IL-1 β Kit
Lot Number: K0082550
Expiration Date: 31 MAR 2026
Catalog Number: K151QP-Series, K156QP-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IL-1 β Antibody	D0082006	2-8°C	31 AUG 2026
Human IL-1 β Plate	Z0048023	2-8°C	28 FEB 2027
Proinflammatory Panel 1 (human) Calibrator Blend	A0080407	2-8°C	31 AUG 2029
Diluent 2	M0020128	\leq -10°C	31 MAR 2026
Diluent 3	M0030129	\leq -10°C	30 NOV 2026

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-1 β	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-1 β	2.14	80% - 120%	< 20%	375	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IL-1 β	546	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	Hamilla, Karen	Quality		30 OCT 2024