

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX and V-PLEX Plus Human IL-1 β Kit**Lot Number:** K0082916**Expiration Date:** 30 SEP 2027**Catalog Number:** K151QP-Series, K156QP-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-1 β Antibody	D0082366	2-8°C	31 MAR 2028
Human IL-1 β 96-Well Small Spot SECTOR Plate	Z0048246	2-8°C	31 AUG 2028
Proinflammatory Panel 1 (human) Calibrator Blend	A0080439	2-8°C	31 JUL 2030
Diluent 2	M0020134	$\leq -10^\circ\text{C}$	30 SEP 2027
Diluent 3	M0030141	$\leq -10^\circ\text{C}$	30 NOV 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-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	592	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		29 APR 2026