

**CERTIFICATE OF ANALYSIS**

**Product Description:** V-PLEX® and V-PLEX Human IL-1 $\beta$  Kit  
**Lot Number:** K0082611  
**Expiration Date:** 31 JAN 2026  
**Catalog Number:** K151QP-Series, K156QP-Series

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IL-1 $\beta$ Antibody	D0081884	2-8°C	31 JAN 2026
Human IL-1 $\beta$ 96-Well Small Spot SECTOR Plate	Z0047973	2-8°C	30 SEP 2026
Proinflammatory Panel 1 (human) Calibrator Blend	A0080378	2-8°C	31 AUG 2028
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 $\beta$	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 $\beta$	2.14	80% - 120%	< 20%	375	80% - 120%	< 20%

**Stock Concentration of Calibrator:**

Spot	Name	Conc.	Units
1	IL-1 $\beta$	627	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
<b>Review/Approval</b>	Karen Hamilla	Quality		06 JAN 2025