

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX PLUS Human IL-1 β Kit
Kit Lot Number: K0081830
Expiration Date: 31 Oct 2023
Kit Catalog Numbers: K151QP-Series; K156QP-Series

Precision and Accuracy 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	600	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-1 β Plate	Z0047583	2-8°C	29 Feb 2024
Anti-hu IL-1 β Antibody	D0081575	2-8°C	31 Aug 2024
Proinflammatory Panel 1 (human) Calibrator Blend	A0080291	2-8°C	30 Apr 2026
Diluent 2	M0020117	\leq -10°C	31 Oct 2023
Diluent 3	M0030109	\leq -10°C	31 Jan 2024

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

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		13 APR 2022