

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IL-1 β Kit
Kit Lot Number: K0081604
Expiration Date: 31 Aug 2022
Kit Catalog Numbers: K151QP-Series, K156QP-Series

Precision and Accuracy Specifications:

		Calibrator		Proinflammatory Panel 1(human) Control 1		Proinflammatory Panel 1(human) Control 2		Proinflammatory Panel 1(human) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-1 β	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

		LLOQ			ULOQ		
Spot	Name	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	605	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-1 β Plate	Z0047429	2-8°C	30 Apr 2023
Anti-hu IL-1 β Antibody	D0081274	2-8°C	31 Aug 2022
Proinflammatory Panel 1 (human) Calibrator Blend	A0080269	2-8°C	30 Sep 2025
Diluent 2	M0020114	$\leq$ -10°C	31 Mar 2023
Diluent 3	M0030104	$\leq$ -10°C	30 Jun 2023

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		11 FEB 2021