

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-1 $\beta$  Kit  
**Kit Lot Number:** K0081561  
**Expiration Date:** 30 Apr 2022  
**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 $\beta$	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 $\beta$	623	pg/ml

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-1 $\beta$ Plate	Z0047401	2-8°C	31 Mar 2023
Anti-hu IL-1 $\beta$ Antibody	D0081182	2-8°C	30 Apr 2022
Proinflammatory Panel 1(human) Calibrator Blend	A0080254	2-8°C	30 Apr 2025
Diluent 2	M0020113	$\leq$ -10°C	30 Nov 2022
Diluent 3	M0030101	$\leq$ -10°C	31 Dec 2022

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	Xiomara Talavera	Quality		16 NOV 2020