

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Mouse IL-1 $\beta$  Kit  
**Kit Lot Number:** K0081756  
**Expiration Date:** 30 Jun 2023  
**Kit Catalog Numbers:** K152QP-Series

**Precision and Accuracy Specifications:**

Spot	Name	Calibrator		Proinflammatory Panel 1 (mouse) Control 1		Proinflammatory Panel 1 (mouse) Control 2		Proinflammatory Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-1 $\beta$	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-1 $\beta$	0.720	75% - 125%	< 25%	1030	75% - 125%	< 25%

**Stock Concentration of Calibrator:**


Spot	Name	Conc.	Units
1	IL-1 $\beta$	1810	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-1 $\beta$ Plate	Z0047548	2-8°C	30 Nov 2023
Anti-ms IL-1 $\beta$ Antibody	D0081439	2-8°C	30 Jun 2023
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080282	2-8°C	31 Mar 2026
Diluent 41	M0410040	$\leq$ -10°C	31 Oct 2023
Diluent 45	M0450039	$\leq$ -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		23 JUL 2021