

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-1 β Kit
Kit Lot Number: K0082396
Expiration Date: 30 Nov 2025
Kit Catalog Numbers: K152QP-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-1 β Plate	Z0047930	2-8°C	30 Apr 2026
Anti-ms IL-1 β Antibody	D0081922	2-8°C	28 Feb 2026
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080351	2-8°C	31 Dec 2027
Diluent 41	M0410053	\leq -10°C	30 Nov 2025
Diluent 45	M0450053	\leq -10°C	31 Mar 2026

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

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 β	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

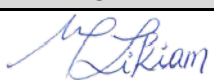
Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-1 β	0.720	75% - 125%	< 25%	1030	75% - 125%	< 25%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IL-1 β	1740	pg/mL

Dilution to Top of Curve = 1X

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

	Name	Function	Signature	Date
Review/Approval	Michelle Likiam	Quality		07 Dec 2023