



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-1β Kit

Lot Number: K0082770

Expiration Date: 30 NOV 2026

Catalog Number: K152QP-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-ms IL-1β Antibody	D0082216	2-8°C	31 JUL 2027
Mouse IL-1β SECTOR Plate	Z0048157	2-8°C	31 JAN 2028
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080401	2-8°C	30 JUN 2029
Diluent 41	M0410058	≤ -10°C	30 SEP 2027
Diluent 45	M0450054	≤ -10°C	30 NOV 2026

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

Precision, Accuracy, and Sensitivity 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-1β	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-1β	0.72	75% - 125%	< 25%	1030	75% - 125%	< 25%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IL-1β	1680	pg/mL

Dilution to top of Curve: 1X

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

	Name	Function	Signature	Date
Review/Approval	Michelle Likiam	Quality		10 OCT 2025