



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-1 β Kit
Lot Number: K0082665
Expiration Date: 30 NOV 2026
Catalog Number: K152QP-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-ms IL-1 β Antibody	D0082216	2-8°C	31 JUL 2027
Mouse IL-1 β 96-Well Small Spot SECTOR Plate	Z0048093	2-8°C	31 AUG 2027
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080373	2-8°C	31 AUG 2028
Diluent 41	M0410056	$\leq -10^{\circ}\text{C}$	31 MAR 2027
Diluent 45	M0450054	$\leq -10^{\circ}\text{C}$	30 NOV 2026

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

Precision, Accuracy, and Sensitivity Specifications:

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

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

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IL-1 β	1700	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	Karen Hamilla	Quality		08 APR 2025