

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-1 β Kit
Kit Lot Number: K0081906
Expiration Date: 31 Jan 2024
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 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	1530	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-1 β Plate	Z0047624	2-8°C	30 Apr 2024
Anti-ms IL-1 β Antibody	D0081534	2-8°C	31 Jan 2024
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080252	2-8°C	30 Apr 2025
Diluent 41	M0410043	\leq -10°C	31 Mar 2024
Diluent 45	M0450043	\leq -10°C	31 Jan 2024

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		06 MAY 2022