



Extension Memo

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IL-1 β Kit
Catalog Number(s): K151QP-Series; K156QP-Series
Lot Number: K0082119
New Expiration Date: 30 Jun 2025

Based on additional testing, the original date of expiration for this kit has been extended by 12 months.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Human IL-1 β Plate	Z0047758	28 Feb 2025	30 Jun 2025
Anti-hu IL-1 β Antibody	D0081575	31 Aug 2024	30 Jun 2025
Diluent 2	M0020120	30 Jun 2024	30 Jun 2025
Diluent 3	M0030113	31 Jul 2024	30 Jun 2025

Function	Name	Signature	Date
Quality	Karen Hamilla		30 Jan 2025

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IL-1 β Kit
Kit Lot Number: K0082119
Expiration Date: 30 Jun 2024
Kit Catalog Numbers: K151QP-Series; K156QP-Series

Precision and Accuracy Specifications:

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

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-1 β	2.14	80% - 120%	< 20%	375	80% - 120%	< 20%

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	IL-1 β	613	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-1 β Plate	Z0047758	2-8°C	28 Feb 2025
Anti-hu IL-1 β Antibody	D0081575	2-8°C	31 Aug 2024
Proinflammatory Panel 1 (human) Calibrator Blend	A0080328	2-8°C	30 Apr 2027
Diluent 2	M0020120	\leq -10°C	30 Jun 2024
Diluent 3	M0030113	\leq -10°C	31 Jul 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		26 SEP 2022