



## Extension Memo

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-12p70 Kit  
**Catalog Number(s):** K151QV-Series  
**Lot Number:** K0082175  
**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-12p70 Plate	Z0047788	31 May 2025	30 Jun 2025
Anti-hu IL-12p70Antibody	D0081643	31 Aug 2024	30 Jun 2025
Diluent 2	M0020120	30 Jun 2024	30 Jun 2025
Diluent 3	M0030114	30 Sep 2024	30 Jun 2025

Function	Name	Signature	Date
Quality	Karen Hamilla		12 Sep 2024

## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-12p70 Kit  
**Kit Lot Number:** K0082175  
**Expiration Date:** 30 Jun 2024  
**Kit Catalog Numbers:** K151QV-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-12p70	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-12p70	1.22	80% - 120%	< 20%	315	80% - 120%	< 20%

**Stock Concentration of Calibrator:**


Spot	Name	Conc.	Units
1	IL-12p70	492	pg/mL

**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-12p70 Plate	Z0047788	2-8°C	31 May 2025
Anti-hu IL-12p70Antibody	D0081643	2-8°C	31 Aug 2024
Proinflammatory Panel 1(human) Calibrator Blend	A0080334	2-8°C	31 Jul 2027
Diluent 2	M0020120	≤ -10°C	30 Jun 2024
Diluent 3	M0030114	≤ -10°C	30 Sep 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		02 JAN 2023