



EXTENSION MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX Plus Human IL-12p70 Kit


Catalog Number(s) / Size(s): K151QV-Series

Lot Number: K0082086

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0082086 has been extended to 30 SEP 2024.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Diluent 2	M0020116	30 Sep 2023	30 SEP 2024
Anti-hu IL-12p70 Antibody	D0081643	31 Aug 2024	30 SEP 2024
Diluent 3	M0030111	31 May 2024	30 SEP 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		19 JUL 2023

This memo supersedes all other statements regarding date of expiration for this product.

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IL-12p70 Kit
Kit Lot Number: K0082086
Expiration Date: 30 Sep 2023
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	473	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-12 p70 Plate	Z0047736	2-8°C	31 Jan 2025
Anti-hu IL-12p70 Antibody	D0081643	2-8°C	31 Aug 2024
Proinflammatory Panel 1 (human) Calibrator Blend	A0080311	2-8°C	31 Oct 2026
Diluent 2	M0020116	≤ -10°C	30 Sep 2023
Diluent 3	M0030111	≤ -10°C	31 May 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		30 SEP 2022