



## Extension Memo

Page 1 of 1

**Product Description:** V-PLEX® and V-PLEX PLUS Human IL-15 Kit  
**Catalog Number(s):** K151RD-Series, K156RD-Series  
**Lot Number:** K0082713  
**New Expiration Date:** 30 Sep 2026

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-15 Plate                         | Z0047851   | 30 Sep 2025              | 30 Sep 2026         |
| Cytokine Panel 1 (human) Calibrator Blend | A0080300   | 31 Jul 2026              | 30 Sep 2026         |
| Anti-hu IL-15 Antibody                    | D0082034   | 31 Aug 2026              | 30 Sep 2026         |

|                 | Name          | Function | Signature                                                                            | Date        |
|-----------------|---------------|----------|--------------------------------------------------------------------------------------|-------------|
| Review/Approval | Karen Hamilla | Quality  |  | 24 Jun 2025 |



## CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX® and V-PLEX PLUS Human IL-15 Kit  
**Lot Number:** K0082713  
**Expiration Date:** 30 SEP 2025  
**Catalog Number:** K151RD-Series; K156RD-Series

**Lot-Specific Components:**

| Description                               | Lot Number | Storage Temperature | Expiration Date |
|-------------------------------------------|------------|---------------------|-----------------|
| Anti-hu IL-15 Antibody                    | D0082034   | 2-8°C               | 31 AUG 2026     |
| Cytokine Panel 1 (human) Calibrator Blend | A0080300   | 2-8°C               | 31 JUL 2026     |
| Human IL-15 Plate                         | Z0047851   | 2-8°C               | 30 SEP 2025     |
| Diluent 3                                 | M0030136   | ≤ -10°C             | 31 MAY 2027     |
| Diluent 43                                | M0430056   | ≤ -10°C             | 30 SEP 2026     |

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

**Precision, Accuracy, and Sensitivity Specifications:**

| Spot | Name  | Calibrator |       | Cytokine Panel 1 (human)<br>Control 1 |       | Cytokine Panel 1 (human)<br>Control 2 |       | Cytokine Panel 1 (human)<br>Control 3 |       |
|------|-------|------------|-------|---------------------------------------|-------|---------------------------------------|-------|---------------------------------------|-------|
|      |       | Recovery   | CV    | Recovery                              | CV    | Recovery                              | CV    | Recovery                              | CV    |
| 1    | IL-15 | 80% - 120% | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% |

| Spot | Name  | LLOQ  |            |       | ULOQ  |            |       |
|------|-------|-------|------------|-------|-------|------------|-------|
|      |       | Value | Recovery   | CV    | Value | Recovery   | CV    |
| 1    | IL-15 | 0.774 | 80% - 120% | < 20% | 525   | 80% - 120% | < 20% |

**Stock Concentration of Calibrator:**

| Spot | Name  | Conc. | Units |
|------|-------|-------|-------|
| 1    | IL-15 | 811   | 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  |  | 23 MAY 2025 |