

**Extension Memo**

**Product Description:** V-PLEX and V-PLEX Plus Cytokine Panel 1 (hu) Kit, QuickPlex  
**Lot Number:** K0082683  
**New Expiration Date:** 30 JUN 2027  
**Catalog Number(s):** K15050-Series, K15057-Series, K151QS-Series, K151RB-Series, K151RC-Series, K151RD-Series, K151RE-Series, K151RF-Series, K151RG-Series, K151RH-Series, K151RI-Series, K151RJ-Series, K156QS-Series, K156RC-Series, K156RD-Series, K156RE-Series, K156RF-Series, K156RG-Series, K156RH-Series, K156RI-Series, K156RJ-Series

Based on additional testing, the original date of expiration for this kit has been extended by 9 months.

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

| Description                  | Lot Number | Original Expiration Date | New Expiration Date |
|------------------------------|------------|--------------------------|---------------------|
| Anti-hu IL-12/23p40 Antibody | D0082137   | 28 FEB 2027              | 30 JUN 2027         |
| Anti-hu IL-5 Antibody        | D0082138   | 28 FEB 2027              | 30 JUN 2027         |
| Diluent 43                   | M0430056   | 30 SEP 2026              | 30 JUN 2027         |
| Diluent 3                    | M0030136   | 31 MAY 2027              | 30 JUN 2027         |

| Function | Name          | Signature                                                                            | Date        |
|----------|---------------|--------------------------------------------------------------------------------------|-------------|
| Quality  | Karen Hamilla |  | 14 Sep 2026 |

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX and V-PLEX Plus Cytokine Panel 1 (hu) Kit, QuickPlex**Lot Number:** K0082683**Expiration Date:** 30 SEP 2026**Catalog Number:** K15050-Series, K15057-Series, K151QS-Series, K151RB-Series, K151RC-Series, K151RD-Series, K151RE-Series, K151RF-Series, K151RG-Series, K151RH-Series, K151RI-Series, K151RJ-Series, K156QS-Series, K156RC-Series, K156RD-Series, K156RE-Series, K156RF-Series, K156RG-Series, K156RH-Series, K156RI-Series, K156RJ-Series**Lot-Specific Components:**

| Description                                                 | Lot Number | Storage Temperature        | Expiration Date |
|-------------------------------------------------------------|------------|----------------------------|-----------------|
| Anti-hu GM-CSF Antibody                                     | D0082279   | 2-8°C                      | 31 DEC 2027     |
| Anti-hu IL-12/23p40 Antibody                                | D0082137   | 2-8°C                      | 28 FEB 2027     |
| Anti-hu IL-15 Antibody                                      | D0082283   | 2-8°C                      | 30 NOV 2027     |
| Anti-hu IL-16 Antibody                                      | D0082182   | 2-8°C                      | 30 JUN 2027     |
| Anti-hu IL-17A Antibody                                     | D0082284   | 2-8°C                      | 30 NOV 2027     |
| Anti-hu IL-1 $\alpha$ Antibody                              | D0082280   | 2-8°C                      | 31 JAN 2028     |
| Anti-hu IL-5 Antibody                                       | D0082138   | 2-8°C                      | 28 FEB 2027     |
| Anti-hu IL-7 Antibody                                       | D0082220   | 2-8°C                      | 31 AUG 2027     |
| Anti-hu TNF- $\beta$ Antibody                               | D0082221   | 2-8°C                      | 31 AUG 2027     |
| Anti-hu VEGF Antibody                                       | D0082439   | 2-8°C                      | 31 AUG 2028     |
| Cytokine Panel 1 (hu) 96 Well 10-Spot QuickPlex Ultra Plate | Z0048063   | 2-8°C                      | 31 JUL 2027     |
| Cytokine Panel 1 (human) Calibrator Blend                   | A0080383   | 2-8°C                      | 31 DEC 2028     |
| Diluent 3                                                   | M0030136   | $\leq -10^{\circ}\text{C}$ | 31 MAY 2027     |
| Diluent 43                                                  | M0430056   | $\leq -10^{\circ}\text{C}$ | 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    | GM-CSF         | 80% - 120% | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% |
| 2    | IL-1 $\alpha$  | 80% - 120% | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% |
| 3    | IL-5           | 80% - 120% | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% |
| 4    | IL-7           | 80% - 120% | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% |
| 5    | IL-12/IL-23p40 | 80% - 120% | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% |
| 6    | IL-15          | 80% - 120% | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% |
| 7    | IL-16          | 80% - 120% | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% |
| 8    | IL-17A         | 80% - 120% | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% |
| 9    | TNF- $\beta$   | 80% - 120% | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% |
| 10   | VEGF-A         | 80% - 120% | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% | 80% - 120%                            | < 20% |



## CERTIFICATE OF ANALYSIS

| Spot | Name           | LLOQ  |            |       | ULOQ  |            |       |
|------|----------------|-------|------------|-------|-------|------------|-------|
|      |                | Value | Recovery   | CV    | Value | Recovery   | CV    |
| 1    | GM-CSF         | 0.842 | 80% - 120% | < 20% | 750   | 80% - 120% | < 20% |
| 2    | IL-1 $\alpha$  | 2.85  | 80% - 120% | < 20% | 278   | 80% - 120% | < 20% |
| 3    | IL-5           | 4.41  | 80% - 120% | < 20% | 562   | 80% - 120% | < 20% |
| 4    | IL-7           | 0.851 | 80% - 120% | < 20% | 563   | 80% - 120% | < 20% |
| 5    | IL-12/IL-23p40 | 1.32  | 80% - 120% | < 20% | 2250  | 80% - 120% | < 20% |
| 6    | IL-15          | 0.774 | 80% - 120% | < 20% | 525   | 80% - 120% | < 20% |
| 7    | IL-16          | 19.1  | 80% - 120% | < 20% | 1870  | 80% - 120% | < 20% |
| 8    | IL-17A         | 3.19  | 80% - 120% | < 20% | 3650  | 80% - 120% | < 20% |
| 9    | TNF- $\beta$   | 0.465 | 80% - 120% | < 20% | 458   | 80% - 120% | < 20% |
| 10   | VEGF-A         | 7.7   | 80% - 120% | < 20% | 562   | 80% - 120% | < 20% |

## Non-Specific Binding Specifications:

| Spot | Name           | Spec   |
|------|----------------|--------|
| 1    | GM-CSF         | < 1.0% |
| 2    | IL-1 $\alpha$  | < 1.0% |
| 3    | IL-5           | < 1.0% |
| 4    | IL-7           | < 1.0% |
| 5    | IL-12/IL-23p40 | < 1.0% |
| 6    | IL-15          | < 1.0% |
| 7    | IL-16          | < 1.0% |
| 8    | IL-17A         | < 1.0% |
| 9    | TNF- $\beta$   | < 1.0% |
| 10   | VEGF-A         | < 1.0% |

## Stock Concentration of Calibrator:

| Spot | Name           | Conc. | Units |
|------|----------------|-------|-------|
| 1    | GM-CSF         | 1110  | pg/mL |
| 2    | IL-1 $\alpha$  | 412   | pg/mL |
| 3    | IL-5           | 826   | pg/mL |
| 4    | IL-7           | 887   | pg/mL |
| 5    | IL-12/IL-23p40 | 3460  | pg/mL |
| 6    | IL-15          | 850   | pg/mL |
| 7    | IL-16          | 2360  | pg/mL |
| 8    | IL-17A         | 6490  | pg/mL |
| 9    | TNF- $\beta$   | 721   | pg/mL |
| 10   | VEGF-A         | 1820  | 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  |  | 03 APR 2025 |