



Extension Memo

Product Description: Diluent 40
Lot Number: M0400022
New Expiration Date: 31 AUG 2027
Catalog Number(s): R50AJ-Series

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

| Function | Name | Signature | Date |
|----------|-----------------|-----------|-------------|
| Quality | Michelle Likiam | | 02 Jun 2026 |



Extension Memo

Product Description: Diluent 42
Lot Number: M0420021
New Expiration Date: 31 AUG 2027
Catalog Number(s): R50AK-Series

Based on additional testing, the original date of expiration for this diluent has been extended by 5 months.

| Function | Name | Signature | Date |
|----------|-----------------|-----------|-------------|
| Quality | Michelle Likiam | | 02 Jun 2026 |

**CERTIFICATE OF ANALYSIS**

Product Description: V-PLEX® and V-PLEX Plus Proinflammatory Panel 2 (rat) Kit
Lot Number: K0082949
Expiration Date: 31 AUG 2027
Catalog Number: K153QP-Series, K153QX-Series, K153QU-Series, K15059-Series, K153QO-Series, K153QR-Series, K153QS-Series, K153QT-Series, K153OD-Series

Lot-Specific Components:

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

| Description | Lot Number | Storage Temperature | Expiration Date |
|--|------------|----------------------------|-----------------|
| Anti-rat IFN- γ Antibody | D0082385 | 2-8°C | 31 JUL 2028 |
| Anti-rat IL-10 Antibody | D0082391 | 2-8°C | 31 JUL 2028 |
| Anti-rat IL-13 Antibody | D0082392 | 2-8°C | 31 MAY 2028 |
| Anti-rat IL-1 β Antibody | D0082387 | 2-8°C | 31 MAY 2028 |
| Anti-rat IL-4 Antibody | D0082386 | 2-8°C | 31 MAY 2028 |
| Anti-rat IL-5 Antibody | D0082388 | 2-8°C | 30 JUN 2028 |
| Anti-rat IL-6 Antibody | D0082389 | 2-8°C | 31 MAY 2028 |
| Anti-rat KC/GRO Antibody | D0082390 | 2-8°C | 30 JUN 2028 |
| Anti-rat TNF- α Antibody | D0082393 | 2-8°C | 30 NOV 2028 |
| Proinflammatory Panel 1 (rat) Calibrator Blend | A0080413 | 2-8°C | 30 APR 2030 |
| Proinflammatory Panel 2 (rat) 96-Well 10-Spot SECTOR Plate | Z0048241 | 2-8°C | 31 AUG 2028 |
| Diluent 40 | M0400022 | $\leq -10^{\circ}\text{C}$ | 31 AUG 2027 |
| Diluent 42 | M0420021 | $\leq -10^{\circ}\text{C}$ | 31 AUG 2027 |
| Blocker H | R0100026 | Room Temperature | 29 FEB 2028 |

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

Precision, Accuracy, and Sensitivity Specifications:

| Spot | Name | Calibrator | | Proinflammatory Panel 1 (rat) Control 1 | | Proinflammatory Panel 1 (rat) Control 2 | | Proinflammatory Panel 1 (rat) Control 3 | |
|------|---------------|------------|-------|---|-------|---|-------|---|-------|
| | | Recovery | CV | Recovery | CV | Recovery | CV | Recovery | CV |
| 1 | IFN- γ | 80% - 120% | < 20% | 75% - 125% | < 25% | 75% - 125% | < 25% | 75% - 125% | < 25% |
| 2 | IL-1 β | 80% - 120% | < 20% | 75% - 125% | < 25% | 75% - 125% | < 25% | 75% - 125% | < 25% |
| 4 | IL-4 | 80% - 120% | < 20% | 75% - 125% | < 25% | 75% - 125% | < 25% | 75% - 125% | < 25% |
| 5 | IL-5 | 80% - 120% | < 20% | 75% - 125% | < 25% | 75% - 125% | < 25% | 75% - 125% | < 25% |
| 6 | IL-6 | 80% - 120% | < 20% | 75% - 125% | < 25% | 75% - 125% | < 25% | 75% - 125% | < 25% |
| 7 | KC/GRO | 80% - 120% | < 20% | 75% - 125% | < 25% | 75% - 125% | < 25% | 75% - 125% | < 25% |
| 8 | IL-10 | 80% - 120% | < 20% | 75% - 125% | < 25% | 75% - 125% | < 25% | 75% - 125% | < 25% |
| 9 | IL-13 | 80% - 120% | < 20% | 75% - 125% | < 25% | 75% - 125% | < 25% | 75% - 125% | < 25% |
| 10 | TNF- α | 80% - 120% | < 20% | 75% - 125% | < 25% | 75% - 125% | < 25% | 75% - 125% | < 25% |



CERTIFICATE OF ANALYSIS

| Spot | Name | LLOQ | | | ULOQ | | |
|------|---------------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | IFN- γ | 39.7 | 75% - 125% | < 25% | 3750 | 75% - 125% | < 25% |
| 2 | IL-1 β | 102 | 75% - 125% | < 25% | 8100 | 75% - 125% | < 25% |
| 4 | IL-4 | 8.00 | 75% - 125% | < 25% | 723 | 75% - 125% | < 25% |
| 5 | IL-5 | 82 | 75% - 125% | < 25% | 3000 | 75% - 125% | < 25% |
| 6 | IL-6 | 96.9 | 75% - 125% | < 25% | 8550 | 75% - 125% | < 25% |
| 7 | KC/GRO | 21.7 | 75% - 125% | < 25% | 728 | 75% - 125% | < 25% |
| 8 | IL-10 | 163 | 75% - 125% | < 25% | 15700 | 75% - 125% | < 25% |
| 9 | IL-13 | 12.5 | 75% - 125% | < 25% | 1080 | 75% - 125% | < 25% |
| 10 | TNF- α | 9.10 | 75% - 125% | < 25% | 793 | 75% - 125% | < 25% |

Non-Specific Binding Specifications:

| Spot | Name | Spec |
|------|---------------|--------|
| 1 | IFN- γ | < 1.0% |
| 2 | IL-1 β | < 1.0% |
| 4 | IL-4 | < 1.0% |
| 5 | IL-5 | < 1.0% |
| 6 | IL-6 | < 1.0% |
| 7 | KC/GRO | < 1.0% |
| 8 | IL-10 | < 1.0% |
| 9 | IL-13 | < 1.0% |
| 10 | TNF- α | < 1.0% |

Stock Concentration of Calibrator:

| Spot | Name | Conc. | Units |
|------|---------------|-------|-------|
| 1 | IFN- γ | 6300 | pg/mL |
| 2 | IL-1 β | 11900 | pg/mL |
| 4 | IL-4 | 1050 | pg/mL |
| 5 | IL-5 | 9540 | pg/mL |
| 6 | IL-6 | 13400 | pg/mL |
| 7 | KC/GRO | 2950 | pg/mL |
| 8 | IL-10 | 23900 | pg/mL |
| 9 | IL-13 | 1670 | pg/mL |
| 10 | TNF- α | 1330 | 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 | | 08 JUN 2026 |