

**CERTIFICATE OF ANALYSIS****Product Description:** V-PLEX® and V-PLEX PLUS Human TARC Gen. B Kit**Lot Number:** K0082635**Expiration Date:** 31 AUG 2026**Catalog Number:** K151ANK-Series; K156ANK-Series**Lot-Specific Components:**

| Description | Lot Number | Storage Temperature | Expiration Date |
|---|------------|---------------------|-----------------|
| Chemokine Panel 1 (human) Gen. B Calibrator | A0080389 | 2-8°C | 30 APR 2029 |
| Human TARC Antibody | D0082258 | 2-8°C | 31 OCT 2027 |
| Human TARC Gen. B Plate | Z0048069 | 2-8°C | 31 JUL 2027 |
| Diluent 3 | M0030128 | ≤ -10°C | 31 OCT 2026 |
| Diluent 57 | M0570015 | ≤ -10°C | 31 AUG 2026 |

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

Precision, Accuracy, and Sensitivity Specifications:

| Spot | Name | Calibrator | | Chemokine Panel 1 (human) Gen. B Control 1 | | Chemokine Panel 1 (human) Gen. B Control 2 | | Chemokine Panel 1 (human) Gen. B Control 3 | |
|------|------------|------------|-------|--|-------|--|-------|--|-------|
| | | Recovery | CV | Recovery | CV | Recovery | CV | Recovery | CV |
| 1 | TARC/CCL17 | 80% - 120% | < 20% | 90% - 130% | < 20% | 90% - 130% | < 20% | 90% - 130% | < 20% |

| Spot | Name | LLOQ | | | ULOQ | | |
|------|------------|-------|------------|-------|-------|------------|-------|
| | | Value | Recovery | CV | Value | Recovery | CV |
| 1 | TARC/CCL17 | 2.14 | 80% - 120% | < 20% | 481 | 75% - 125% | < 20% |

Stock Concentration of Calibrator:

| Spot | Name | Conc. | Units |
|------|------------|-------|-------|
| 1 | TARC/CCL17 | 778 | 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 | | 17 MAR 2025 |