

**CERTIFICATE OF ANALYSIS**

**Product Description:** V-PLEX® and V-PLEX PLUS Human TARC Gen. B Kit  
**Lot Number:** K0082352  
**Expiration Date:** 31 JUL 2025  
**Catalog Number:** K151ANK-Series; K156ANK-Series

**Lot-Specific Components:**

| Description                                       | Lot Number | Storage Temperature | Expiration Date |
|---|------------|---------------------|-----------------|
| Chemokine Panel 1 (human) Gen. B Calibrator Blend | A0080377   | 2-8°C               | 31 AUG 2028     |
| Human TARC Antibody                               | D0082132   | 2-8°C               | 31 JAN 2027     |
| Human TARC Gen. B Plate                           | Z0047903   | 2-8°C               | 28 FEB 2026     |
| Diluent 3   | M0030123   | ≤ -10°C             | 31 JAN 2026     |
| Diluent 57  | M0570011   | ≤ -10°C             | 31 JUL 2025     |

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% | 80% - 120%                                 | < 20% | 80% - 120%                                 | < 20% | 80% - 120%                                 | < 20% |

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

**Stock Concentration of Calibrator:**

| Spot | Name       | Conc. | Units |
|------|------------|-------|-------|
| 1    | TARC/CCL17 | 744   | 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        |
|------------------------|---------------|----------|-----------|-------------|
| <b>Review/Approval</b> | Karen Hamilla | Quality  |           | 30 MAY 2024 |