



## CERTIFICATE OF ANALYSIS

**Product Description:** V-PLEX and V-PLEX Plex Plus Ab40 Peptide (4G8) Kit**Lot Number:** K0082738**Expiration Date:** 30 APR 2027**Catalog Number:** K150SJ-Series**Lot-Specific Components:**

| Description                                      | Lot Number | Storage Temperature        | Expiration Date |
|--------------------------------------------------|------------|----------------------------|-----------------|
| Anti-A $\beta$ 4G8 Antibody                      | D0082419   | 2-8°C                      | 31 JUL 2028     |
| A $\beta$ 40 Peptide 96-Well 4-Spot SECTOR Plate | Z0057592   | 2-8°C                      | 30 NOV 2027     |
| A $\beta$ 40 Blocker                             | R0010925   | 2-8°C                      | 30 SEP 2027     |
| Diluent 100                                      | Y0270344   | 2-8°C                      | 30 APR 2027     |
| Diluent 35                                       | M0350039   | 2-8°C                      | 30 JUN 2027     |
| A $\beta$ 1-40 Peptide (4G8)                     | A0080425   | $\leq -70^{\circ}\text{C}$ | 31 AUG 2027     |

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

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

|      |              | Calibrator |       | Neurodegeneration Control 1 (4G8) |       | Neurodegeneration Control 2 (4G8) |       | Neurodegeneration Control 3 (4G8) |       |
|------|--------------|------------|-------|-----------------------------------|-------|-----------------------------------|-------|-----------------------------------|-------|
| Spot | Name         | Recovery   | CV    | Recovery                          | CV    | Recovery                          | CV    | Recovery                          | CV    |
| 1    | A $\beta$ 40 | 80% - 120% | < 20% | 75% - 125%                        | < 20% | 75% - 125%                        | < 20% | 75% - 125%                        | < 20% |

|      |              | LLOQ  |            |       | ULOQ  |            |       |
|------|--------------|-------|------------|-------|-------|------------|-------|
| Spot | Name         | Value | Recovery   | CV    | Value | Recovery   | CV    |
| 1    | A $\beta$ 40 | 20    | 80% - 120% | < 20% | 6000  | 80% - 120% | < 20% |

**Stock Concentration of Calibrator:**

| Spot | Name         | Conc.  | Units |
|------|--------------|--------|-------|
| 1    | A $\beta$ 40 | 616000 | pg/mL |

Dilution to top of Curve: 40X

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

|                 | Name          | Function | Signature                                                                            | Date        |
|-----------------|---------------|----------|--------------------------------------------------------------------------------------|-------------|
| Review/Approval | Karen Hamilla | Quality  |  | 18 AUG 2025 |