



## CERTIFICATE OF ANALYSIS

**Product Description:** U-PLEX® and U-PLEX Plus Human  $\alpha$ -Synuclein Kit  
**Lot Number:** K0082509  
**Expiration Date:** 31 JUL 2026  
**Catalog Number:** K151WK-Series

**Lot-Specific Components:**

| Description  | Lot Number | Storage Temperature | Expiration Date |
|--|------------|---------------------|-----------------|
| MSD GOLD™ 96-well Small Spot Streptavidin SECTOR Plate | Z0022421   | 2-8°C               | 31 DEC 2026     |
| Diluent 49   | M0490019   | ≤ -10°C             | 31 JAN 2027     |
| Anti-hu $\alpha$ -Synuclein                            | D00U2357   | ≤ -70°C             | 31 DEC 2026     |
| Biotin Anti-hu $\alpha$ -Synuclein                     | B00U2401   | ≤ -70°C             | 31 JUL 2026     |
| Human $\alpha$ -Synuclein Calibrator                   | A0080399   | ≤ -70°C             | 30 NOV 2026     |

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

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

| Spot | Name                | Calibrator |       | Human $\alpha$ -Synuclein Control 1 |       | Human $\alpha$ -Synuclein Control 2 |       | Human $\alpha$ -Synuclein Control 3 |       |
|------|---------------------|------------|-------|-------------------------------------|-------|-------------------------------------|-------|-------------------------------------|-------|
|      |                     | Recovery   | CV    | Recovery                            | CV    | Recovery                            | CV    | Recovery                            | CV    |
| 1    | $\alpha$ -Synuclein | 80% - 120% | < 20% | 80% - 120%                          | < 20% | 80% - 120%                          | < 20% | 80% - 120%                          | < 20% |

| Spot | Name                | LLOQ  |            |       | ULOQ  |            |       |
|------|---------------------|-------|------------|-------|-------|------------|-------|
|      |                     | Value | Recovery   | CV    | Value | Recovery   | CV    |
| 1    | $\alpha$ -Synuclein | 8     | 80% - 120% | < 20% | 6800  | 80% - 120% | < 20% |

**Stock Concentration of Calibrator:**

| Spot | Name                | Conc.  | Units |
|------|---------------------|--------|-------|
| 1    | $\alpha$ -Synuclein | 210000 | pg/mL |

Dilution to top of Curve: 20X

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  |           | 07 OCT 2024 |