



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: A β 1-42 Peptide (6E10)

Catalog Number: C01LB-2

Lot Number: A0080200

Expiration Date: 30JUN2021

Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration of Materials Tested:

| | Assay |
|-----------------------|-----------------------|
| | A β 1-42 (6E10) |
| Calibrator (pg/mL) | 70 000 |

Concentration: 40X

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

| | Name | Function | Signature | Date |
|------------------------|--------------|-------------------|---------------------|-------------|
| Review/Approval | Dawn Stollar | Quality Assurance | <i>Dawn Stollar</i> | 28 JAN 2019 |