



## CERTIFICATE OF ANALYSIS


**Product Description:** A $\beta$ 1-42 Peptide (4G8)  
**Catalog Number:** C003A-2  
**Lot Number:** A0080426  
**Expiration Date:** 31 AUG 2027  
**Storage Temperature:**  $\leq -70^{\circ}\text{C}$

**Concentration Assignment:**

Analyte	Stock Concentration	Unit of Measure
A $\beta$ 42	79200	pg/mL

**Dilution to Top of Curve:** 40 X

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. 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		03 APR 2025