



CERTIFICATE OF ANALYSIS

Product Description: A β 1-42 Peptide (4G8)**Catalog Number:** C003A-2**Lot Number:** A0080471**Expiration Date:** 31 JAN 2029**Storage Temperature:** $\leq -70^{\circ}\text{C}$ **Concentration Assignment:**

Analyte	Stock Concentration	Unit of Measure
A β 42	82400	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
Review/Approval	Karen Hamilla	Quality		31 AUG 2026