



CERTIFICATE OF ANALYSIS

Product Description: A β 1-40 Peptide (4G8)
Catalog Number: C002Z-2
Lot Number: A0080449
Expiration Date: 31 AUG 2028
Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
A β 40(4G8)	576000	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		20 MAR 2026