



## CERTIFICATE OF ANALYSIS

**Product Description:** A $\beta$ 1-40 Peptide (4G8)  
**Catalog Number:** C002Z-2  
**Lot Number:** A0080359  
**Expiration Date:** 31 JUL 2025  
**Storage Temperature:**  $\leq -70^{\circ}\text{C}$

## Concentration Assignment

Analyte	Stock Concentration	Unit of Measure
A $\beta$ 1-40	616000	pg/mL

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

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		24 MAR 2023