



## CERTIFICATE OF ANALYSIS


**Product Description:** A $\beta$ 1-38 Peptide (6E10)  
**Catalog Number:** C00NZ-2  
**Lot Number:** A0080451  
**Expiration Date:** 30 APR 2028  
**Storage Temperature:**  $\leq -70^{\circ}\text{C}$

**Concentration Assignment:**

Analyte	Stock Concentration	Unit of Measure
A $\beta$ 38	552000	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		21 NOV 2025