



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** A $\beta$ 1-38 Peptide (6E10)

**Catalog Number:** C00NZ-2

**Lot Number:** A0080287

**Expiration Date:** 30 SEP 2023


**Storage Temperature:**  $\leq -70^{\circ}\text{C}$

### Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
A $\beta$ 38	488000	pg/mL

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

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		28 SEP 2021