



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: A β 1-38 Peptide (4G8)

Catalog Number: C002Y-2

Lot Number: A0080335

Expiration Date: 31 JAN 2025

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

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
A β 38	544000	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	<i>Karen Hamilla</i>	08 SEP 2022