



**CERTIFICATE OF ANALYSIS**

Page 1 of 1

**Product Description:** A $\beta$ 1-38 Peptide (4G8)

**Catalog Number:** C002Y-2

**Lot Number:** A0080292

**Expiration Date:** 31 OCT 2023

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

**Concentration Assignment:**

Analyte	Stock Concentration	Unit of Measure
A $\beta$ 1-38	472000	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		02 NOV 2021