



**RFE MEMO**


Page 1 of 1

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

**Catalog Number(s) / Size(s):** C002Y-2

**Lot Number:** A0080189

Based on additional testing, the original date of expiration for this calibrator has been extended by 12 months. The expiration date of A0080189 has been extended to 30 SEP 2022.

Function	Name	Signature	Date
Quality	Karen Hamilla		10 JUN 2021

This memo supersedes all other statements regarding date of expiration for this product.



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C002Y-2

**Lot Number:** A0080189

**Expiration Date:** 30 SEP 2021

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

### Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
A $\beta$ 1-38 (4G8)	480 000	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	Dawn Stollar	Quality		30 APR 2019