



RFE MEMO


Page 1 of 1

Product Description(s): A β 1-38 Peptide

Catalog Number: C00NZ-2

Lot Number: A0080181

Based on additional testing, the original date of expiration for this Calibrator has been extended by 12 months. The expiration date of A0080181 has been extended to 28 FEB 2022.

Function	Name	Signature	Date
Quality	Karen Hamilla		04 MAY 2021

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



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: A β 1-38 Peptide

Catalog Number: C00NZ-2

Lot Number: A0080181

Expiration Date: 28 FEB 2021

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

Concentration of Materials Tested:

	Assay
	A β 1-38
Calibrator (pg/mL)	516 000

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	09 OCT 2018