



RFE MEMO


Page 1 of 1

Product Description(s): A β 1-42 Peptide (4G8)

Catalog Number(s) / Size(s): C003A-2

Lot Number: A0080201

Based on additional testing, the original date of expiration for this calibrator has been extended by 10 months. The expiration date of A0080201 has been extended to 31 MAY 2022.

Function	Name	Signature	Date
Quality	Karen Hamilla		08 APR 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 1-42 Peptide (4G8)

Catalog Number: C003A-2

Lot Number: A0080201

Expiration Date: 31 JUL 2021


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

Concentration of Materials Tested:

	Assay
	A β 1-42 Peptide (4G8)
Calibrator (pg/mL)	99 200

Concentration: 40X

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		15 APR 2019