



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: A β 1-40 Peptide

Catalog Number: C000A-2

Lot Number: A0080190

Expiration Date: 30 APR 2021

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

Concentration of Materials Tested:

	Assay
	A β 1-40
Calibrator (pg/mL)	604 000

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>	11 DEC 2018