



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: A β 1-42 Peptide

Catalog Number: C01LB-2

Lot Number: A0080140

Expiration Date: 30 NOV 2019

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

Concentration of Materials Tested:

	Assay
	A β 1-42 Peptide
Calibrator (pg/mL)	94 400

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>	15 AUG 2017