



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: A β 1-38 Peptide

Catalog Number: C00NZ-2

Lot Number: A0080134

Expiration Date: 30 NOV 2019

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

Concentration of Materials Tested:

	Assay
	A β 1-38
Calibrator (pg/mL)	564 000

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

	Name	Function	Signature	Date
Review/Approval	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	27JUL2017