



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07049

Product Description: A β 1-38 Peptide
Catalog Number: C00NZ-2
Lot Number: A0040165
Expiration Date: 30 Sep 2018
Storage Temperature: $\leq -70^{\circ}\text{C}$

Stock Concentration of Materials Tested:

	Assay
	A β 38
Calibrator (pg/mL)	339 200

Statement: The above product was manufactured and tested using MSD procedures and it conforms to MSD's specifications. 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>	02 AUG 2016