



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** A $\beta$ 1-38 Peptide

**Catalog Number:** C00NZ-2

**Lot Number:** A0080148

**Expiration Date:** 31 AUG 2020

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

### Concentration of Materials Tested:

	Assay
	A $\beta$ 1-38
Calibrator (pg/mL)	592 000

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	24 APR 2018