



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: A β 1-42 Peptide (6E10)

Catalog Number: C01LB-2

Lot Number: A0080245

Expiration Date: 31JUL 2022


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

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
A β 1-42 (6E10)	73600	pg/ml

Dilution to Top of Curve: 40X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		21 FEB 2020