



CERTIFICATE OF ANALYSIS

Product Description: Aβ1-42 Peptide (6E10)
Catalog Number: C01LB-2
Lot Number: A0080409
Expiration Date: 28 FEB 2027
Storage Temperature: ≤ -70°C

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
Aβ42	72800	pg/mL

Dilution to Top of Curve: 40 X

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality	<i>Karen Hamilla</i>	02 OCT 2024