



CERTIFICATE OF ANALYSIS

Control Pack Description: Neurodegeneration Control Pack 1 (6E10)
Catalog Number: C41LB-1
Lot Number: CP000187
Expiration Date: 31 DEC 2028

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Neurodegeneration Control 1 (6E10)	A00C1041	≤ -70°C	31 DEC 2028
Neurodegeneration Control 2 (6E10)	A00C1042	≤ -70°C	31 DEC 2028
Neurodegeneration Control 3 (6E10)	A00C1043	≤ -70°C	31 DEC 2028

Concentration Assignment:

Analyte	Control 1	Control 2	Control 3	Unit of Measure
Aβ40	9220	3090	524	pg/mL
Aβ38	4130	991	218	pg/mL
Aβ42	1450	473	187	pg/mL

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		18 FEB 2026