

ALZpath Simoa® p-Tau 217 v2 EQC Kit Certificate of Testing



Data below represents results generated on Simoa Analyzers.

Lot Number:	999025
Item #:	104372
Manufacturing Date:	06-Dec-2023
Expiration Date:	N/A
Platform(s):	HD-X

Control Materials**				** Ranges shown here are for reference only. We recommend customer to generate their own lot-specific control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
EQC 1	999013	0.556	0.389 - 0.723	
EQC 2	999014	0.284	0.199 - 0.369	

Review/Approval

Director, Accelerator Services

Title

L. Ni 12 Dec 2023

Signature/Date