

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



Data below represents results generated on Simoa Analyzers.

Lot Number:	999062
Item #:	104372
Manufacturing Date:	14-Mar-2024
Expiration Date:	N/A
Platform(s):	HD-X

Control Materials**			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
EQC 1	999041	1.31	0.955 - 1.77
EQC 2	999042	0.247	0.189 - 0.350

** Ranges shown here are for reference only. We recommend customer to generate their own lot-specific control ranges.

Review/Approval

Director, Accelerator Services

Title

L. J. 26 Mar 2024

Signature/Date