

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



Data below represents results generated on Simoa Analyzers.

Lot Number:	999009
Item #:	104372
Manufacturing Date:	10-Jul-2023
Expiration Date:	N/A
Platform(s):	HD-X

Control Materials**			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
EQC 1	999006	2.15	1.51 - 2.80
EQC 2	999007	0.242	0.169 - 0.314

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

Review/Approval

Director, Accelerator Services

Title

[Handwritten Signature] 28 Jul 2023

Signature/Date