

# ALZpath Simoa® p-Tau 217 v2 Assay Kit Certificate of Analysis

Lot Number:	999115
Item #:	104371
Manufacturing Date:	18-Mar-2025
Expiration Date:	18-Mar-2026
Platform(s):	HD-X



Data below represents results generated on Simoa Analyzers.

Component	Select one				Lot Number
	Reagent	Stock	Conc.	Ref. Cal.	
Bead	<input checked="" type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	999111
Detector	<input checked="" type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	999112
SβG	<input checked="" type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	999113
Calibrator	<input type="checkbox"/>	<input type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>	999109
Sample Diluent	N/A				999114
Calibrator Diluent	N/A				999114
Bead Diluent	N/A				N/A
Detector Diluent	N/A				N/A
SβG Diluent	N/A				N/A
RGP Reagent (Part # 103159)	N/A				N/A

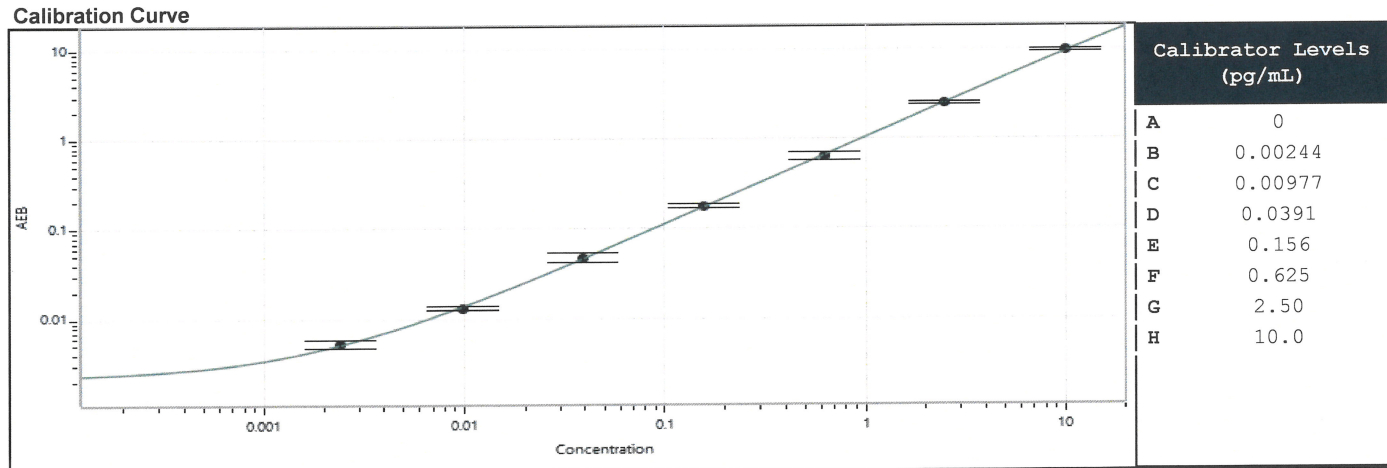
\* Unless specified, RGP is not Kit Lot Specific

Control Materials**			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	N/A	N/A	N/A
Control 2	N/A	N/A	N/A

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

Assigned value for Calibrator Stock Concentration (pg/mL)	1,060	Not Applicable	<input type="checkbox"/>
LOD (Limit of Detection). See TECH-0163.		Not Provided	<input checked="" type="checkbox"/>
LLOQ (Lower Limit of Quantitation). See TECH-0163.		Not Provided	<input checked="" type="checkbox"/>

Calibration Curve



Review/Approval

Director, Accelerator Services  
Title

*[Signature]* 26 Mar 2025  
Signature/Date