

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

Lot Number:	999024
Item #:	104371
Manufacturing Date:	26-Sep-2023
Expiration Date:	N/A
Platform(s):	HD-X



Data below represents results generated on Simoa Analyzers.

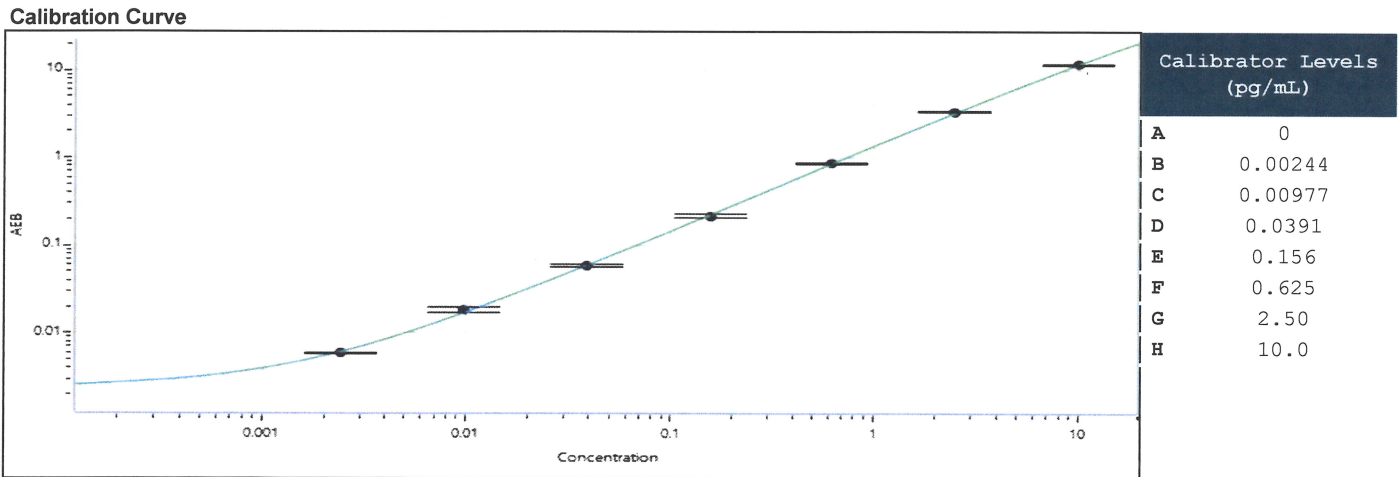
Component	Select one		Lot Number
Bead	Reagent	<input checked="" type="checkbox"/> Stock	999016
Detector	Reagent	<input checked="" type="checkbox"/> Stock	999017
SβG	Reagent	<input checked="" type="checkbox"/> Stock	999018
Calibrator	Ref. Cal.	<input type="checkbox"/> Conc. <input checked="" type="checkbox"/>	999012
Sample Diluent	N/A		999019
Calibrator Diluent	N/A		999019
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**				** 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)	
Control 1	N/A	N/A	N/A	
Control 2	N/A	N/A	N/A	

Assigned value for Calibrator Stock Concentration (pg/mL)	917	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] 29 Sep 2023

Signature/Date