

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

Lot Number:	999036
Item #:	104371
Manufacturing Date:	06-Dec-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	999032
Detector	Reagent	<input checked="" type="checkbox"/> Stock	999033
SβG	Reagent	<input checked="" type="checkbox"/> Stock	999034
Calibrator	Ref. Cal.	<input type="checkbox"/> Conc. <input checked="" type="checkbox"/>	999028
Sample Diluent	N/A		999035
Calibrator Diluent	N/A		999035
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

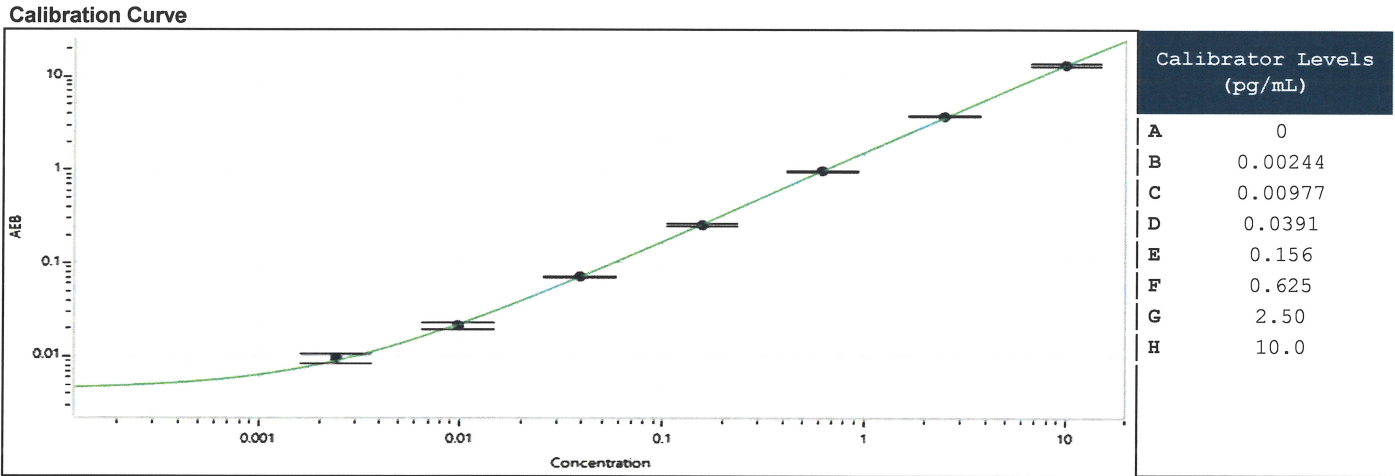
* Unless specified, RGP is not Kit Lot Specific

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)	1,110	Not Applicable	
LOD (Limit of Detection). See TECH-0163.		Not Provided	x
LLOQ (Lower Limit of Quantitation). See TECH-0163.		Not Provided	x

Calibration Curve



Review/Approval

Director, Accelerator Services

Title

Signature/Date

[Handwritten Signature] 12 Dec 2023