

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

Lot Number:	999008
Item #:	104371
Manufacturing Date:	10-Jul-2023
Expiration Date:	N/A
Platform(s):	HD-X



Data below represents results generated on Simoa Analyzers.

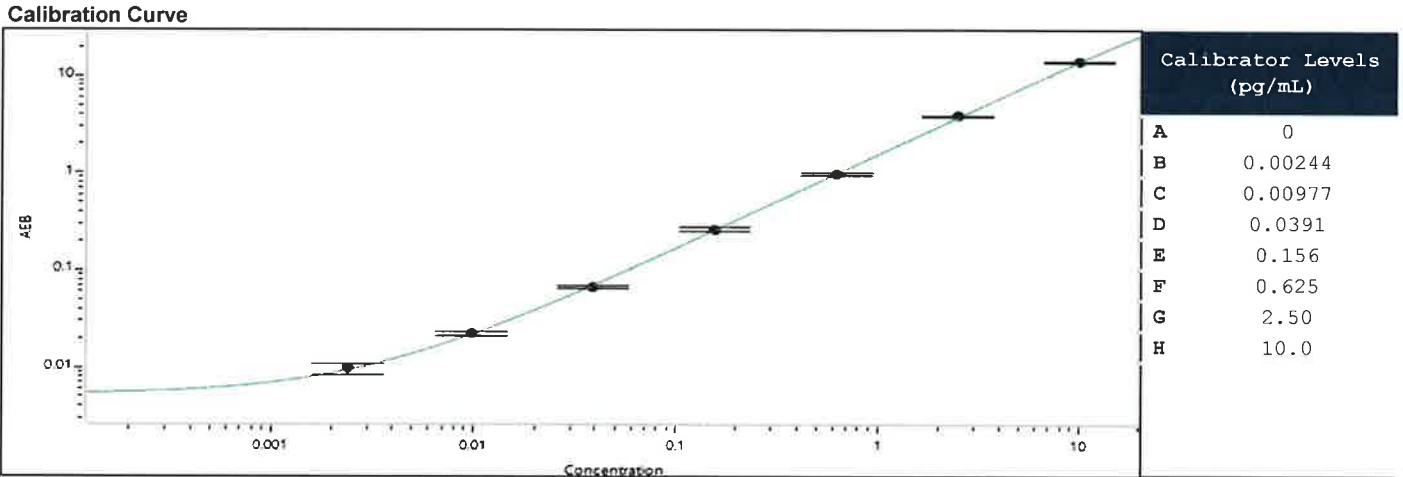
Component	Select one			Lot Number	
Bead	Reagent	<input checked="" type="checkbox"/>	Stock	999001	
Detector	Reagent	<input checked="" type="checkbox"/>	Stock	999002	
SPG	Reagent	<input checked="" type="checkbox"/>	Stock	999003	
Calibrator	Ref. Cal.	<input type="checkbox"/>	Conc.	<input checked="" type="checkbox"/>	999005
Sample Diluent	N/A			999004	
Calibrator Diluent	N/A			999004	
Bead Diluent	N/A			N/A	
Detector Diluent	N/A			N/A	
SPG 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)	335	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/Date

[Handwritten Signature] 7/28/2023