

ALZpath Simoa® p-Tau 217 v2 Assay Kit

Lot Number:	MAB231122
Item #:	104371
Manufacturing Date:	30-Nov-22
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 <input type="checkbox"/>	MAB061222B
Detector	Reagent <input checked="" type="checkbox"/> Stock <input type="checkbox"/>	MAB061222D
SβG	Reagent <input checked="" type="checkbox"/> Stock <input type="checkbox"/>	MAB061222E
Calibrator	Ref. Cal <input checked="" type="checkbox"/> Conc. <input type="checkbox"/>	MAB221122
Sample Diluent	N/A	MAB061222F
Calibrator Diluent	N/A	MAB061222F
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

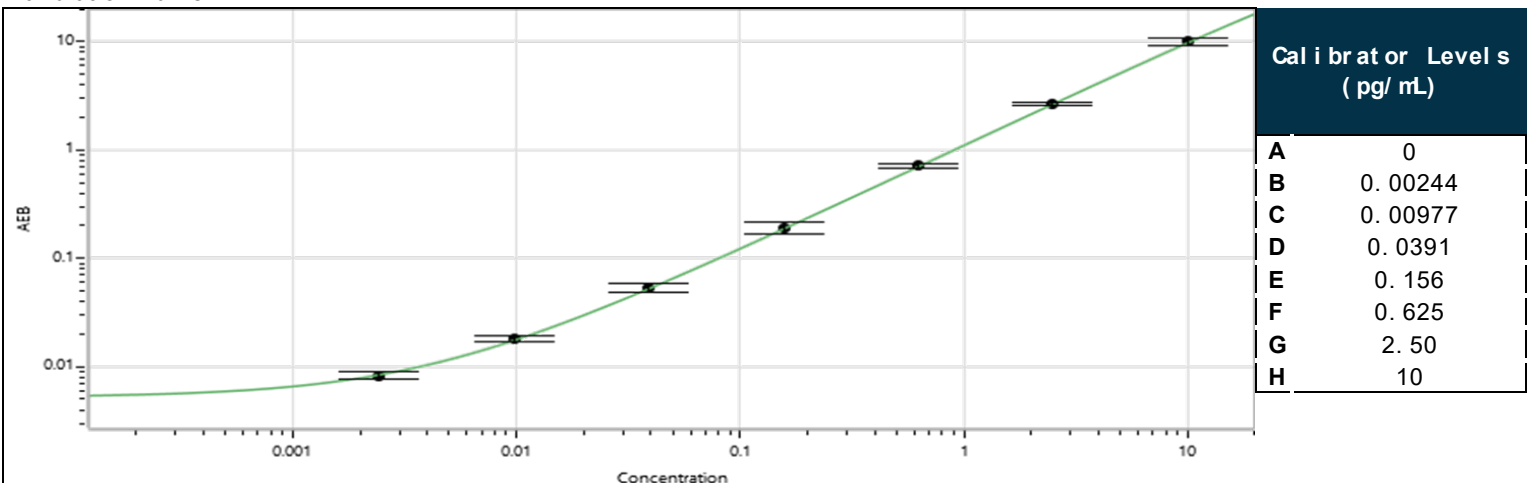
* Unless specified, RGP is not Kit Lot Specific

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)	1,130	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

Hong Ni 16 Jun 2023

Title

Signature/Date