

Product Number:	104112
Lot Number:	504236
Paired Reagent Kit: (104111)	504132
Expiration:	12-Aug-2025
Platform(s):	SR-X and HD-1/HD-X



Component	Part Number	Lot Number
pTau-181 V2.1 Sample Diluent Reagent	104101	408112
pTau-181 V2.1 Calibrator Diluent Reagent	104103	404322
pTau-181 V2.1 Calibrator Concentrate	104102	411405

Data below represents results generated on the Simoa® HD-X Analyzer.

Calibrator Concentrate Value Assignment (pg/mL)	36311
--	-------

Review/Approval

Cody Carlson

QC Manager

[Handwritten Signature] 25 JUN 2024

Name

Title

Signature/Date