

Product Number:	104798
Lot Number:	504402
Paired Reagent Kit Lot Number:	504346
Expiration:	4-Apr-2029
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Calibrator Concentrate	104793	409518
Control 1 Concentrate	104795	409519
Control 2 Concentrate	104796	409520

Calibrator Concentrate Reconstitution Volume (uL)	1751
Control 1 Concentrate Reconstitution Volume (uL)	1751
Control 2 Concentrate Reconstitution Volume (uL)	1751

Review/Approval

TAM LEE DOC CONTROL II, QA Manager 23 OCT 2024
 Name Title Signature/Date