

Product Number:	104650
Lot Number:	504458
Paired Reagent Kit Lot Number:	504238
Expiration:	9-Jan-2029
Platform(s):	HD-X

Component	Part Number	Lot Number
Calibrator Concentrate	104614	400907
Control 1 Concentrate	104616	400908
Control 2 Concentrate	104617	400909



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

Calibrator Concentrate Reconstitution Volume (uL)	1517
Control 1 Concentrate Reconstitution Volume (uL)	841
Control 2 Concentrate Reconstitution Volume (uL)	841

Review/Approval

TAM LEE DOC CONTROL II, QA Tam Lee 20 NOV 2024
 Name Title Signature/Date