

Product Number:	104571
Lot Number:	504565
Paired Reagent Kit Lot Number:	504500
Expiration:	21-Feb-2029
Platform(s):	HD-X and SR-X



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

Component	Part Number	Lot Number
Calibrator Concentrate	104566	405210
Control 1 Concentrate	104568	405211
Control 2 Concentrate	104569	405212

Calibrator Concentrate Reconstitution Volume (uL)	943
Control 1 Concentrate Reconstitution Volume (uL)	943
Control 2 Concentrate Reconstitution Volume (uL)	943

Review/Approval

Kyle Giroux Senior Document Control Specialist, QA [Signature] 27Mar2025
 Name Title Signature/Date