

Product Number:	104571
Lot Number:	536621
Paired Reagent Kit Lot Number:	504600
Expiration:	21-Feb-2029
Platform(s):	HD-X and SR-X

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



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

Calibrator Concentrate Reconstitution Volume (uL)	766
Control 1 Concentrate Reconstitution Volume (uL)	766
Control 2 Concentrate Reconstitution Volume (uL)	766

Review/Approval

Dhvani Patel QA Specialist II [Signature] 10 Nov 2025
 Name Title Signature/Date