

Product Number:	105797
Lot Number:	504695
Paired Reagent Kit Lot Number:	504628
Expiration:	20-Mar-2030
Platform(s):	HD-X

Component	Part Number	Lot Number
Calibrator Concentrate	105792	507910
Control 1 Concentrate	105794	507911
Control 2 Concentrate	105795	507912



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

Calibrator Concentrate Reconstitution Volume (uL)	1303
Control 1 Concentrate Reconstitution Volume (uL)	500
Control 2 Concentrate Reconstitution Volume (uL)	500

Review/Approval

Dhwani Patel QA Specialist II, QA DP | 19 AUG 2025
 Name Title Signature/Date