

Product Number:	105738
Lot Number:	504667
Paired Reagent Kit Lot Number:	504577
Expiration:	13-Feb-2030
Platform(s):	HD-X

Component	Part Number	Lot Number	Control Lot Number
Calibrator Concentrate	105733	504410	504410A
Control 1 Concentrate	105735	504412	504412A
Control 2 Concentrate	105736	504413	504413A



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

Calibrator Concentrate Reconstitution Volume (uL)	762
Control 1 Concentrate Reconstitution Volume (uL)	762
Control 2 Concentrate Reconstitution Volume (uL)	300

Review/Approval

D. Kesh Patel Sr. Doc. Control - QA ET 11/24/2025
Name Title Signature/Date