

Product Number:	104543
Lot Number:	504743
Paired Reagent Kit Lot Number:	504497
Expiration:	9-Jan-2029
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number
Calibrator Concentrate	104538	400911
Control 1 Concentrate	104540	400912
Control 2 Concentrate	104541	400913



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

Calibrator Concentrate Reconstitution Volume (uL)	1177
Control 1 Concentrate Reconstitution Volume (uL)	975
Control 2 Concentrate Reconstitution Volume (uL)	975

Review/Approval

Dhruvi Patel QA Specialist II, QA DP / 01 Oct 2025
 Name Title Signature/Date