

Product Number:	105635
Lot Number:	504668
Paired Reagent Kit Lot Number:	504580
Expiration:	31-Oct-2029
Platform(s):	HD-X

Component	Part Number	Lot Number
Calibrator Concentrate	105630	430514
Control 1 Concentrate	105632	430516
Control 2 Concentrate	105633	430517



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

Calibrator Concentrate Reconstitution Volume (uL)	676
Control 1 Concentrate Reconstitution Volume (uL)	676
Control 2 Concentrate Reconstitution Volume (uL)	676

Review/Approval

Dilish Patel
Name

Sr. Doc. Control - QA
Title

AT 16/11/2025
Signature/Date