

Product Number:	104543
Lot Number:	504370
Paired Reagent Kit Lot Number:	504279
Expiration:	5-Dec-2028
Platform(s):	HD-X

Component	Part Number	Lot Number
Calibrator Concentrate	104538	333907
Control 1 Concentrate	104540	333909
Control 2 Concentrate	104541	333910



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

Calibrator Concentrate Reconstitution Volume (uL)	2000
Control 1 Concentrate Reconstitution Volume (uL)	2000
Control 2 Concentrate Reconstitution Volume (uL)	2000

Review/Approval

Dipesh Patel
Name

Doc control II-QA
Title

[Signature] / 10 Sep 2024
Signature/Date