

Product Number:	104628
Lot Number:	504347
Paired Reagent Kit Lot Number:	503994
Expiration:	21-Jun-2028
Platform(s):	HD-X

Component	Part Number	Lot Number
Calibrator Concentrate	104361	317309
Control 1 Concentrate	104362	317310
Control 2 Concentrate	104363	317311



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

Calibrator Concentrate Reconstitution Volume (uL)	1236
Control 1 Concentrate Reconstitution Volume (uL)	1236
Control 2 Concentrate Reconstitution Volume (uL)	1236

Review/Approval

Robert Somogyi

QC Manager

[Handwritten Signature] 14 AUG 2024

Name

Title

Signature/Date