

Product Number:	103837
Lot Number:	504367
Paired Reagent Kit Lot Number:	504314
Expiration:	30-Oct-2029
Platform(s):	HD-X

Component	Part Number	Lot Number
Calibrator Concentrate	104443	330312
Control 1 Concentrate	104445	330314
Control 2 Concentrate	104446	330315



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

Calibrator Concentrate Reconstitution Volume (uL)	914
Control 1 Concentrate Reconstitution Volume (uL)	914
Control 2 Concentrate Reconstitution Volume (uL)	914

Review/Approval

*Dhvani Patel*

*DOC Control I QA*

*DPP / 12 Sep 2024*

Name

Title

Signature/Date