

Product Number:	104589
Lot Number:	504454
Paired Reagent Kit Lot Number:	504407
Expiration:	23-Oct-2029
Platform(s):	HD-X

Component	Part Number	Lot Number
Calibrator Concentrate	104584	329610
Control 1 Concentrate	104586	329612
Control 2 Concentrate	104587	329613



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

Calibrator Concentrate Reconstitution Volume (uL)	806
Control 1 Concentrate Reconstitution Volume (uL)	806
Control 2 Concentrate Reconstitution Volume (uL)	806

Review/Approval

Shwami Patel
Name

DOC Control I, QA
Title

SP 14 Nov 2024
Signature/Date