

Product Number:	104956
Lot Number:	504671
Paired Reagent Kit Lot Number:	504427
Expiration:	10-Jun-2029
Platform(s):	HD-X

Component	Part Number	Lot Number
Calibrator Concentrate	104951	416213
Control 1 Concentrate	104953	416215
Control 2 Concentrate	104954	416216



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

Calibrator Concentrate Reconstitution Volume (uL)	940
Control 1 Concentrate Reconstitution Volume (uL)	940
Control 2 Concentrate Reconstitution Volume (uL)	940

Review/Approval

Dhruvi Patel

Name

Doc Control II, QA

Title

SDP / 17 Jul 2025

Signature/Date