

Product Number:	104798
Lot Number:	504740
Paired Reagent Kit Lot Number:	504275
Expiration:	4-Apr-2029
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number
Calibrator Concentrate	104793	409512
Control 1 Concentrate	104795	409513
Control 2 Concentrate	104796	409514



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

Calibrator Concentrate Reconstitution Volume (uL)	1465
Control 1 Concentrate Reconstitution Volume (uL)	1465
Control 2 Concentrate Reconstitution Volume (uL)	1465

Review/Approval

Dhruvi Patel QA Specialist II, QA DP | 30 Sep 2025

Name Title Signature/Date