

Product Number:	104692
Lot Number:	504585
Paired Reagent Kit Lot Number:	504379
Expiration:	29-Jan-2029
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number
Calibrator Concentrate	104687	402944
Control 1 Concentrate	104689	402945
Control 2 Concentrate	104690	402946

Calibrator Concentrate Reconstitution Volume (uL)	1037
Control 1 Concentrate Reconstitution Volume (uL)	1037
Control 2 Concentrate Reconstitution Volume (uL)	1037



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

Review/Approval

Dhvani Patel Doc Control / II, QA  24 Jul 2025
 Name Title Signature/Date