

Product Number:	104692
Lot Number:	536642
Paired Reagent Kit Lot Number:	504630
Expiration:	8-Feb-2029
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number
Calibrator Concentrate	104687	403918
Control 1 Concentrate	104689	403919
Control 2 Concentrate	104690	403920



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

Calibrator Concentrate Reconstitution Volume (uL)	1393
Control 1 Concentrate Reconstitution Volume (uL)	1393
Control 2 Concentrate Reconstitution Volume (uL)	1393

Review/Approval

Dharmi Patel QA Specialist II DP / 01 Dec 2025
 Name Title Signature/Date