

Product Number:	104429
Lot Number:	504368
Paired Reagent Kit Lot Number:	504269
Expiration:	2-Nov-2029
Platform(s):	HD-X

Component	Part Number	Lot Number
Calibrator Concentrate	104483	330609
Control 1 Concentrate	104426	330611
Control 2 Concentrate	104427	330612



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

Calibrator Concentrate Reconstitution Volume (uL)	756
Control 1 Concentrate Reconstitution Volume (uL)	756
Control 2 Concentrate Reconstitution Volume (uL)	756

Review/Approval

Dhruvini Patel / Doc Control / I QA
 Name Title

DPP / 11 Sep 2024
 Signature/Date