

Product Number:	104628
Lot Number:	504437
Paired Reagent Kit Lot Number:	504231
Expiration:	21-Jun-2028
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Calibrator Concentrate	104361	404519
Control 1 Concentrate	104362	404520
Control 2 Concentrate	104363	404521

Calibrator Concentrate Reconstitution Volume (uL)	1480
Control 1 Concentrate Reconstitution Volume (uL)	1480
Control 2 Concentrate Reconstitution Volume (uL)	1480

Review/Approval

TAM LEE DOC CONTROL #, QA Tam Lee 23 OCT 2024

Name

Title

Signature/Date