

Product Number:	104692
Lot Number:	504436
Paired Reagent Kit Lot Number:	504342
Expiration:	29-Jan-2029
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Calibrator Concentrate	104687	402909
Control 1 Concentrate	104689	402910
Control 2 Concentrate	104690	402911

Calibrator Concentrate Reconstitution Volume (uL)	911
Control 1 Concentrate Reconstitution Volume (uL)	911
Control 2 Concentrate Reconstitution Volume (uL)	911

Review/Approval

TAM LEE DOC CONTROLL II, QA Tam Lee 23 OCT 2024  
 Name Title Signature/Date