

Product Number:	105067
Lot Number:	504430
Paired Reagent Kit Lot Number:	504364
Expiration:	23-Apr-2029
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Calibrator Concentrate	105062	412220
Control 1 Concentrate	105064	412221
Control 2 Concentrate	105065	412222

Calibrator Concentrate Reconstitution Volume (uL)	1000
Control 1 Concentrate Reconstitution Volume (uL)	1000
Control 2 Concentrate Reconstitution Volume (uL)	1000

Review/Approval

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