

Product Number:	104671
Lot Number:	504464
Paired Reagent Kit Lot Number:	504383
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	104666	402948
Control 1 Concentrate	104668	402949
Control 2 Concentrate	104669	402950

Calibrator Concentrate Reconstitution Volume (uL)	1144
Control 1 Concentrate Reconstitution Volume (uL)	1144
Control 2 Concentrate Reconstitution Volume (uL)	1144

Review/Approval

TAM LEE DOC CONTROL, QA Nam Lee 05 DEC 2024
 Name Title Signature/Date