

Product Number:	104823
Lot Number:	504373
Paired Reagent Kit Lot Number:	504356
Expiration:	22-Jun-2028
Platform(s):	HD-X



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

Component	Part Number	Lot Number	Control Lot Number
Calibrator Concentrate	104818	317324	317324A
Control 1 Concentrate	104820	317325	317325A
Control 2 Concentrate	104821	317326	317326A

Calibrator Concentrate Reconstitution Volume (uL)	1350
Control 1 Concentrate Reconstitution Volume (uL)	1350
Control 2 Concentrate Reconstitution Volume (uL)	1350

Review/Approval

TAM LEE DOC CONTROL II, QA      Tam Lee 20 NOV 2024

Name

Title

Signature/Date