

Product Number:	104544
Lot Number:	504366
Paired Reagent Kit Lot Number:	504182
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	104466	402917
Control 1 Concentrate	104468	402918
Control 2 Concentrate	104469	402919

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

Review/Approval

*Dhruvi Patel*

*Doc Control I, QA*

*EDP 03 Oct 2024*

Name

Title

Signature/Date