

Product Number:	104513
Lot Number:	504369
Paired Reagent Kit Lot Number:	504084
Expiration:	2-Nov-2029
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Calibrator Concentrate	104508	330613
Control 1 Concentrate	104510	330615
Control 2 Concentrate	104511	330616

Calibrator Concentrate Reconstitution Volume (uL)	788
Control 1 Concentrate Reconstitution Volume (uL)	788
Control 2 Concentrate Reconstitution Volume (uL)	788

Review/Approval

*Dhvani Patel*

*Doc Control I QA*

*DPP 13 Sep 2024*

Name

Title

Signature/Date