

Product Number:	104513
Lot Number:	536633
Paired Reagent Kit Lot Number:	504321
Expiration:	2-Nov-2029
Platform(s):	HD-X and SR-X

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



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

Calibrator Concentrate Reconstitution Volume (uL)	811
Control 1 Concentrate Reconstitution Volume (uL)	811
Control 2 Concentrate Reconstitution Volume (uL)	811

Review/Approval

Dhwani Patel QA specialist II *DP* 02 Dec 2025

Name

Title

Signature/Date