

Product Number:	104692
Lot Number:	504585
Paired Reagent Kit Lot Number:	504476
Expiration:	29-Jan-2029
Platform(s):	HD-X

Component	Part Number	Lot Number
Calibrator Concentrate	104687	402944
Control 1 Concentrate	104689	402945
Control 2 Concentrate	104690	402946



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

Calibrator Concentrate Reconstitution Volume (uL)	1053
Control 1 Concentrate Reconstitution Volume (uL)	1053
Control 2 Concentrate Reconstitution Volume (uL)	1053

Review/Approval

William O Quality Systems Engineer - I

Name

Title

William O 03APR2025

Signature/Date