

Product Number:	104571
Lot Number:	504637
Paired Reagent Kit Lot Number:	504536
Expiration:	21-Feb-2029
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number
Calibrator Concentrate	104566	405210
Control 1 Concentrate	104568	405211
Control 2 Concentrate	104569	405212



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

Calibrator Concentrate Reconstitution Volume (uL)	871
Control 1 Concentrate Reconstitution Volume (uL)	871
Control 2 Concentrate Reconstitution Volume (uL)	871

Review/Approval

Dipesh Patel

Name

Sr. Dir. Control-QA

Title

24 / 11 Jun 2025

Signature/Date