

Product Number:	104571
Lot Number:	504710
Paired Reagent Kit Lot Number:	504576
Expiration:	21-Feb-2029
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number
Calibrator Concentrate	104566	406408
Control 1 Concentrate	104568	406409
Control 2 Concentrate	104569	406410



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

Calibrator Concentrate Reconstitution Volume (uL)	868
Control 1 Concentrate Reconstitution Volume (uL)	868
Control 2 Concentrate Reconstitution Volume (uL)	868

Review/Approval

Dhwani Patel

QA Specialist II, QA

CDP | *03 Sep 2025*

Name

Title

Signature/Date