

Product Number:	104571
Lot Number:	504377
Paired Reagent Kit Lot Number:	504180
Expiration:	11-Jul-2028
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Calibrator Concentrate	104566	406702
Control 1 Concentrate	104568	406703
Control 2 Concentrate	104569	406704

Calibrator Concentrate Reconstitution Volume (uL)	649
Control 1 Concentrate Reconstitution Volume (uL)	649
Control 2 Concentrate Reconstitution Volume (uL)	649

Review/Approval

Dhruvi Patel

DOC Control I QA

DPP 13 Sep 2024

Name

Title

Signature/Date