

Product Number:	104628
Lot Number:	504530
Paired Reagent Kit Lot Number:	504501
Expiration:	21-Jun-2028
Platform(s):	HD-X and SR-X



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

Component	Part Number	Lot Number
Calibrator Concentrate	104361	404519
Control 1 Concentrate	104362	404520
Control 2 Concentrate	104363	404521

Calibrator Concentrate Reconstitution Volume (uL)	1387
Control 1 Concentrate Reconstitution Volume (uL)	1387
Control 2 Concentrate Reconstitution Volume (uL)	1387

Review/Approval

Dhwami Patel

DOC Control I, QA

SDP 18 Feb 2025

Name

Title

Signature/Date