

Product Number:	104628
Lot Number:	504535
Paired Reagent Kit Lot Number:	504534
Expiration:	21-Jun-2028
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number
Calibrator Concentrate	104361	412713
Control 1 Concentrate	104362	412714
Control 2 Concentrate	104363	412715



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

Calibrator Concentrate Reconstitution Volume (uL)	1388
Control 1 Concentrate Reconstitution Volume (uL)	1388
Control 2 Concentrate Reconstitution Volume (uL)	1388

Review/Approval

Dipesh Patel

Name

Ssr. Doc. Control-QA

Title

CT 01/14/2025

Signature/Date