

Product Number:	104544
Lot Number:	536645
Paired Reagent Kit Lot Number:	504676
Expiration:	21-Jun-2028
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number
Calibrator Concentrate	104466	317331
Control 1 Concentrate	104468	317355
Control 2 Concentrate	104469	317356



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

Calibrator Concentrate Reconstitution Volume (uL)	972
Control 1 Concentrate Reconstitution Volume (uL)	972
Control 2 Concentrate Reconstitution Volume (uL)	972

Review/Approval

Dhruvi Patel

QA Specialist II

DP / 01 Dec 2025

Name

Title

Signature/Date