

Product Number:	104671
Lot Number:	536654
Paired Reagent Kit Lot Number:	504629
Expiration:	22-Jun-2028
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number	Control Lot Number
Calibrator Concentrate	104666	317341	N/A
Control 1 Concentrate	104668	317342	N/A
Control 2 Concentrate	104669	317343	317343A



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

Calibrator Concentrate Reconstitution Volume (uL)	1144
Control 1 Concentrate Reconstitution Volume (uL)	1144
Control 2 Concentrate Reconstitution Volume (uL)	1144

Review/Approval

William Chan

document control specialist I

William Chan 04DEC2025

Name

Title

Signature/Date