

Product Number:	104628
Lot Number:	504573
Paired Reagent Kit Lot Number:	504511
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)	1201
Control 1 Concentrate Reconstitution Volume (uL)	1201
Control 2 Concentrate Reconstitution Volume (uL)	1201

Review/Approval

William O
Quality Systems Engineer I
William O / 02 APR 2025

Name
Title
Signature/Date