

<b>Product Number:</b>	104628
<b>Lot Number:</b>	504316
<b>Paired Reagent Kit Lot Number:</b>	504147
<b>Expiration:</b>	21-Jun-2028
<b>Platform(s):</b>	HD-X

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



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

<b>Calibrator Concentrate Reconstitution Volume (uL)</b>	1387
<b>Control 1 Concentrate Reconstitution Volume (uL)</b>	1387
<b>Control 2 Concentrate Reconstitution Volume (uL)</b>	1387

Review/Approval

Robert Somogyi

QC Manager

*[Handwritten Signature]* 17 JUL 2024

Name

Title

Signature/Date