

<b>Product Number:</b>	104905
<b>Lot Number:</b>	504463
<b>Paired Reagent Kit Lot Number:</b>	504411
<b>Expiration:</b>	25-Apr-2029
<b>Platform(s):</b>	HD-X



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

Component	Part Number	Lot Number
Calibrator Concentrate	104900	411616
Control 1 Concentrate	104902	411618
Control 2 Concentrate	104903	411619

Calibrator Concentrate Reconstitution Volume (uL)	488
Control 1 Concentrate Reconstitution Volume (uL)	488
Control 2 Concentrate Reconstitution Volume (uL)	488

Review/Approval

TAM LEE      DOC CONTROL II, QA      Tam Lee 05 DEC 2024  
 Name                      Title                      Signature/Date