

Product Number:	104570
Lot Number:	504735
Expiration:	2-Jan-2027
Platform(s):	HD-X and SR-X

Component	Part Number	Lot Number
Bead Reagent	104546	518407
Detector Reagent	104548	519001
SBG Reagent	104550	521905
Sample Diluent	104552	518308
Calibrator Diluent Reagent	104745	518309
Control Diluent Reagent	104554	518310
Calibrator Concentrate	104566	405210
Control 1 Concentrate	104568	405211
Control 2 Concentrate	104569	405212
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific



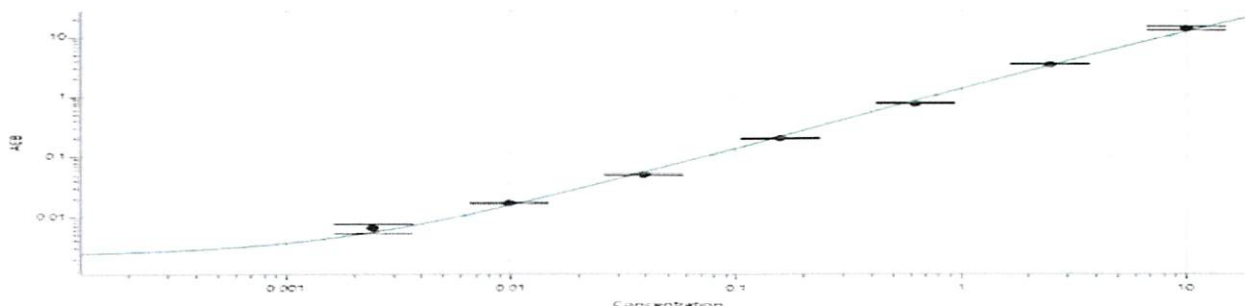
Data below represents results generated on the Simoa® HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	406409	0.758	0.517 - 0.961
Control 2 Concentrate	406410	8.06	5.15 - 9.56

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Reconstitution Volume (uL)	729
Control 1 Concentrate Reconstitution Volume (uL)	729
Control 2 Concentrate Reconstitution Volume (uL)	729

Calibration Curve



Calibrator Levels (pg/mL)

A	0
B	0.00244
C	0.00977
D	0.0391
E	0.156
F	0.625
G	2.50
H	10.0

Review/Approval

Dharmi Patel QA Specialist II, QA *SDR/30SEP2025*

Name

Title

Signature/Date

September 26, 2024

Subject: Simoa® Advantage PLUS Control Concentrate Lots

Dear Valued Customer,

We would like to inform you about a minor detail in the packaging of your Advantage PLUS kit. The kit you received contains Control Concentrates (Level 1 and Level 2) with a different lot number than the one used internally for lot release testing. Both lot numbers refer to materials from the same master batch, and the difference simply reflects different shipment dates. There is no impact on the quality or performance of the product.

If you have any further questions or concerns, please don't hesitate to contact Quanterix's technical support team at TechSupport@Quanterix.com.

Sincerely,

A handwritten signature in black ink, appearing to read 'B. Ciccariello', with a long horizontal stroke extending to the right.

Brian Ciccariello | VP, Quality & Regulatory Affairs