

Product Number:	104428
Lot Number:	504357
Expiration:	19-Feb-2026
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	104409	423208
Detector Reagent	104411	423304
SBG Reagent	104413	423305
Sample Diluent	104415	423222
Calibrator Diluent	104418	423220
Control Diluent	104476	423221
Calibrator Concentrate	104483	412912
Control 1 Concentrate	104426	412910
Control 2 Concentrate	104427	412911
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

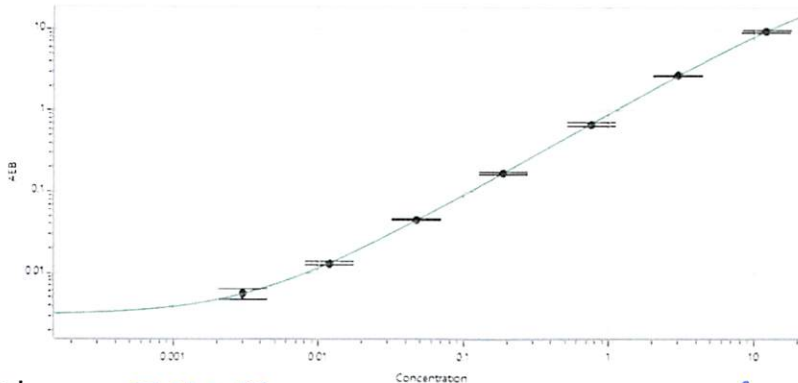
¹ Reagents are not Kit Lot Specific

Release Materials*			* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	
Control 1 Concentrate	330611	0.204	0.117 - 0.236
Control 2 Concentrate	330612	5.94	3.89 - 7.22

Calibrator Concentrate Reconstitution Volume (uL)	832
Control 1 Concentrate Reconstitution Volume (uL)	832
Control 2 Concentrate Reconstitution Volume (uL)	832

Calibrator Levels (pg/mL)	
A	0.000
B	0.003
C	0.012
D	0.047
E	0.188
F	0.750
G	3.00
H	12.0

Calibration Curve



Review/Approval

TAM LEE DOC CONTROL 11, QA Tam Lee 20 SEP 2024
 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 blue ink, appearing to read 'B. Ciccariello', with a long horizontal stroke extending to the right.

Brian Ciccariello | VP, Quality & Regulatory Affairs