

Product Number:	104797
Lot Number:	504346
Expiration:	12-Feb-2026
Platform(s):	HD-X



Component	Part Number	Lot Number
Bead Reagent	104778	422506
Detector Reagent	104780	422601
SBG Reagent	104782	422602
Plasma Sample Diluent	104783	422517
CSF Sample Diluent	104784	422520
Calibrator Diluent	104786	422617
Control Diluent	104787	422618
Calibrator Concentrate	104793	409518
Control 1 Concentrate	104795	409519
Control 2 Concentrate	104796	409520
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

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

¹ Reagents are not Kit Lot Specific

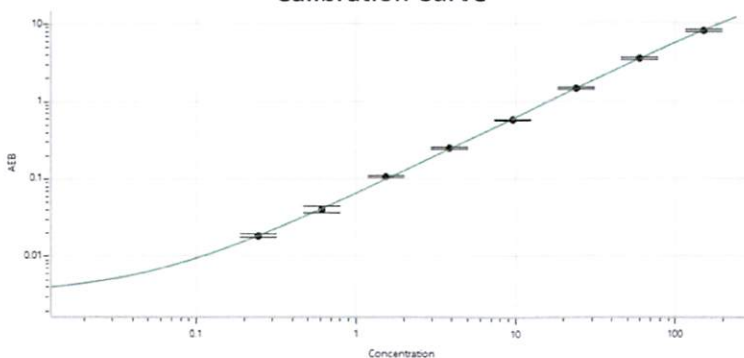
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	409513	7.96	5.25 - 9.87
Control 2 Concentrate	409514	293	212 - 394

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

Calibrator Concentrate Reconstitution Volume (uL)	1751
Control 1 Concentrate Reconstitution Volume (uL)	1751
Control 2 Concentrate Reconstitution Volume (uL)	1751

Calibrator Levels (pg/mL)	
A	0.000
B	0.246
C	0.614
D	1.54
E	3.84
F	9.60
G	24.0
H	60.0
I	150

Calibration Curve



Review/Approval

Dipesh Patel *Doc control II QA*
Name Title

[Signature] *02 OCT 2024*
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".

Brian Ciccariello | VP, Quality & Regulatory Affairs