

Product Number:	103836
Lot Number:	504314
Expiration:	11-Jan-2026
Platform(s):	HD-X



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

Component	Part Number	Lot Number
Bead Reagent	103820	419702
Detector Reagent	103822	419804
SBG Reagent	103824	419805
Sample Diluent	103826	419316
Calibrator Diluent	103829	419314
Control Diluent	104435	419315
Calibrator Concentrate	104443	330312
Control 1 Concentrate	104445	330314
Control 2 Concentrate	104446	330315
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific

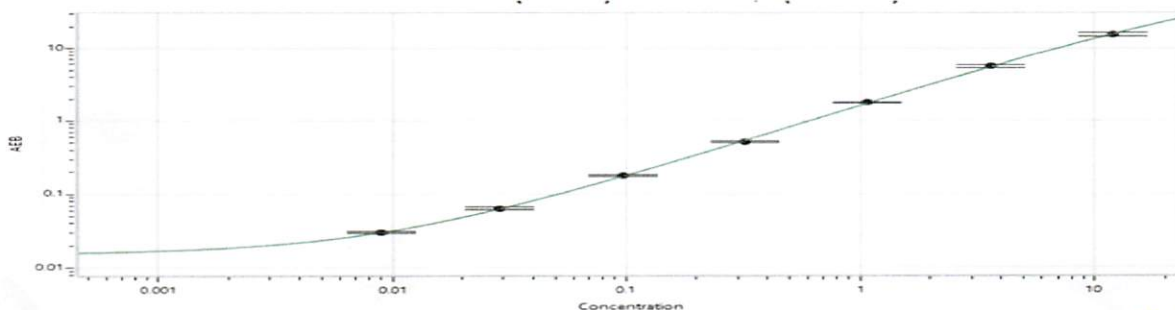
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	330314	0.194	0.135 - 0.251
Control 2 Concentrate	330315	3.53	2.74 - 5.08

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

Calibrator Concentrate Reconstitution Volume (uL)	914
Control 1 Concentrate Reconstitution Volume (uL)	914
Control 2 Concentrate Reconstitution Volume (uL)	914

Calibrator Levels (pg/mL)	
A	0
B	0.009
C	0.029
D	0.098
E	0.325
F	1.08
G	3.60
H	12.0

Calibration Curve



Review/Approval

Cody Carlson

QC Manager

Name

Title

[Signature] 09AUG 2024
Signature/Date