

Product Number:	103277
Lot Number:	504308
Expiration:	5-Sep-2025
Platform(s):	HD-1/HD-X



Component	Part Number	Lot Number
Bead Reagent	103265	418703
Detector Reagent	103268	418704
SBG Reagent	103269	418705
Sample Diluent	103270	418706
Calibrator Diluent	103266	418710
Calibrator Concentrate	103272	420401
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

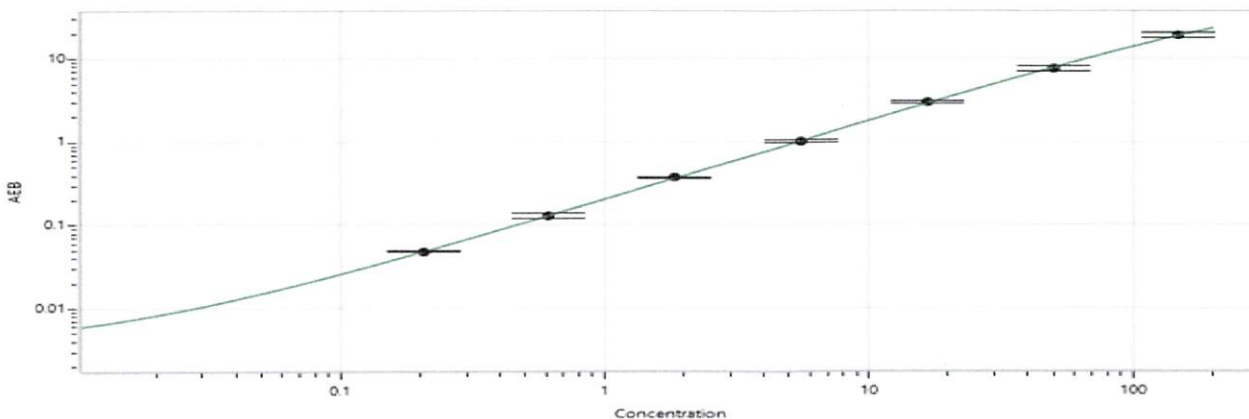
Data below represents results generated on the Simoa® HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	420401	3.58	3.00 - 5.00
Control 2	420401	96.5	75.0 - 125

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

Calibrator Concentrate Value Assignment (pg/mL)	4933
--	------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.206
C	0.617
D	1.85
E	5.56
F	16.7
G	50.0
H	150

Review/Approval

Cody Carlson

QC Manager

Name

Title

[Signature] 06 AUG 2024
Signature/Date