

Product Number:	101622
Lot Number:	504247
Expiration:	14-Feb-2025
Platform(s):	SR-X and HD-1/HD-X



Component	Part Number	Lot Number
Bead Reagent	104233	324408
Detector Reagent	104347	325101
SBG Reagent	104349	325004
Sample Diluent	104351	323617
Calibrator Diluent	101627	322623
Calibrator Concentrate	100652	325803
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

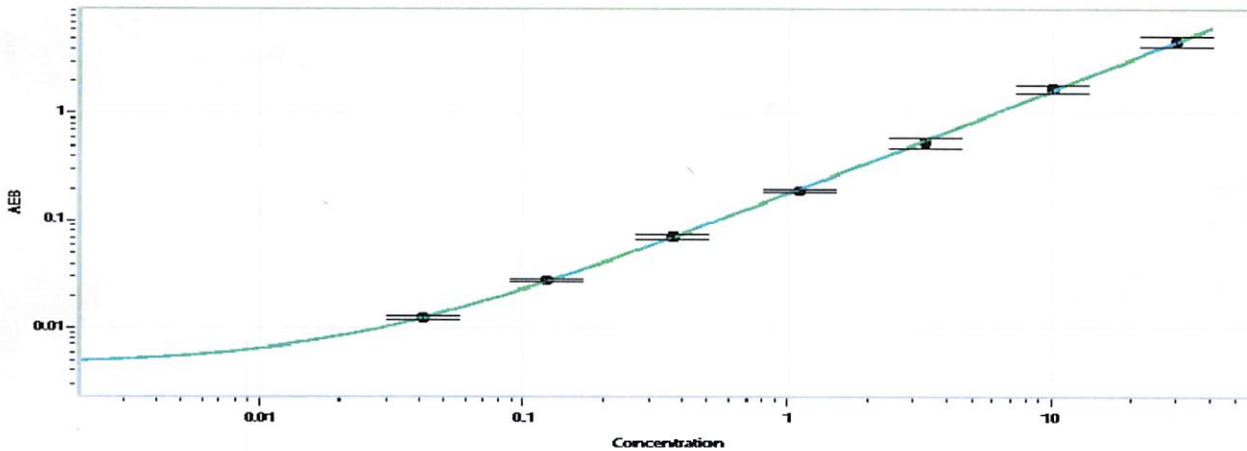
Data below represents results generated on the Simoa® HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	108406	1.56	1.23 - 1.92
Control 2	108406	21.7	17.5 - 28.9
Panel 1	029301	55.8	39.8 - 59.7
Panel 2	029301	6.17	4.35 - 6.53
Panel 3	029301	1.50	1.22 - 1.82

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

Calibrator Concentrate Value Assignment (pg/mL)	492
--	-----

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Cody Carlson
Name Title
QC Manager

[Signature] 26 APR 2024
Signature/Date