

Product Number:	102215
Lot Number:	504130
Expiration:	19-Jul-2025
Platform(s):	SR-X and HD-1/HD-X



Component	Part Number	Lot Number
Bead Reagent	104297	403605
Detector Reagent	104338	403114
SBG Reagent	104340	403606
Sample Diluent	104342	403607
Calibrators	104343	403601
Control 1	104344	403602
Control 2	104345	403602
RGP Reagent	103159	N/A ¹

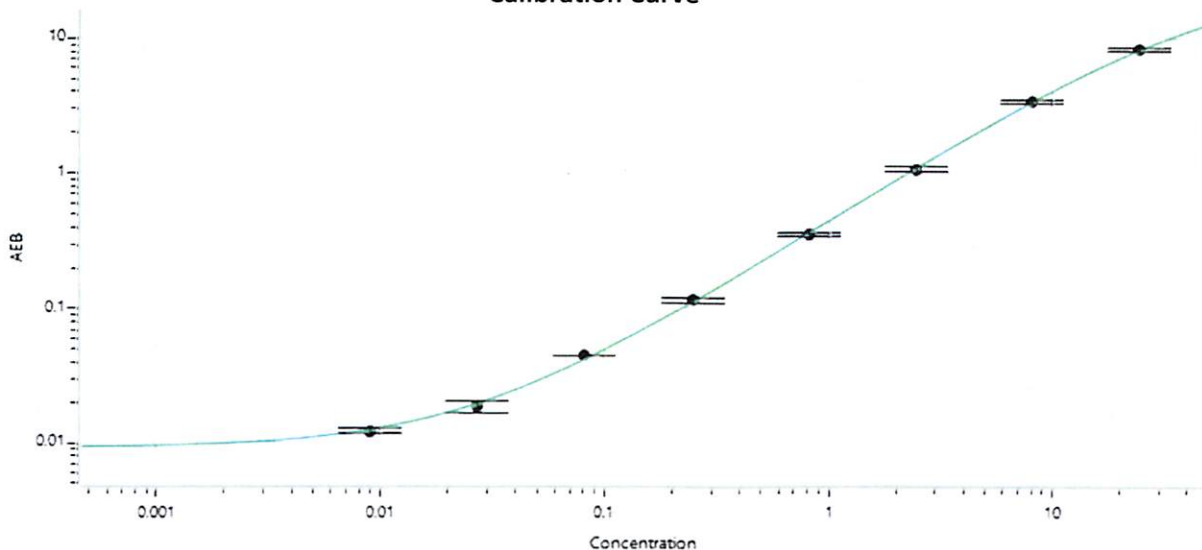
¹ Reagents are not Kit Lot Specific

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

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	403602	0.084	0.057 - 0.100
Control 2	403602	7.86	5.84 - 9.73
Panel 1	019601	0.188	0.127 - 0.243
Panel 2	019601	1.94	1.56 - 2.34
Panel 3	019601	11.0	9.44 - 14.2

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.009
C	0.027
D	0.081
E	0.250
F	0.811
G	2.45
H	8.09
I	24.4

Review/Approval

Cody Carlson

QC Manager

[Signature] 28 FEB 2024

Name

Title

Signature/Date