

<b>Product Number:</b>	104111
<b>Lot Number:</b>	504045
<b>Expiration Date:</b>	18-Jul-25
<b>Platform(s):</b>	SR-X and HD-1/HD-X



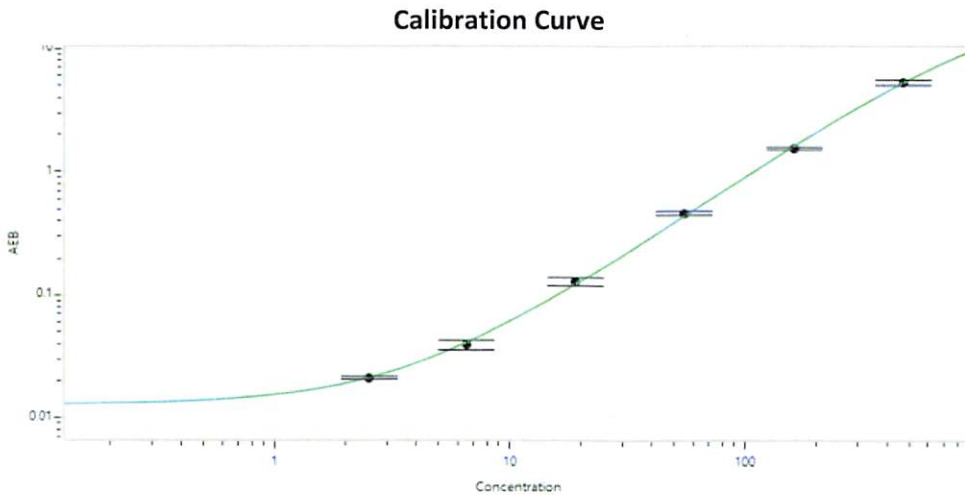
Component	Part Number	Lot Number
Bead Reagent	104098	401809
Detector Reagent	104099	401903
SBG Reagent	104100	401904
Sample Diluent	104101	401908
Calibrators	104106	404301
Control 1	104107	404302
Control 2	104108	404302
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> 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	404302	46.4	36.3 - 54.5
Control 2	404302	1156	919 - 1379
Panel 1	301702	17.0	12.7 - 20.9
Panel 2	301702	14.5	12.3 - 20.9
Panel 3	309704	465	331 - 645
Panel 4	406103	207	158 - 236

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



Calibrator Levels (pg/mL)	
A	0.000
B	2.53
C	6.53
D	19.0
E	55.2
F	162
G	464

Review/Approval

Cody Carlson

QC Manager

*Cody Carlson* 20MAR2024  
Signature/Date

Name

Title