

Product Number:	104111
Lot Number:	504313
Expiration Date:	14-Apr-25
Platform(s):	SR-X and HD-1/HD-X



Component	Part Number	Lot Number
Bead Reagent	104098	328705
Detector Reagent	104099	329108
SBG Reagent	104100	329109
Sample Diluent	104101	329110
Calibrators	104106	333803
Control 1	104107	333804
Control 2	104108	333804
RGP Reagent	103159	N/A ¹

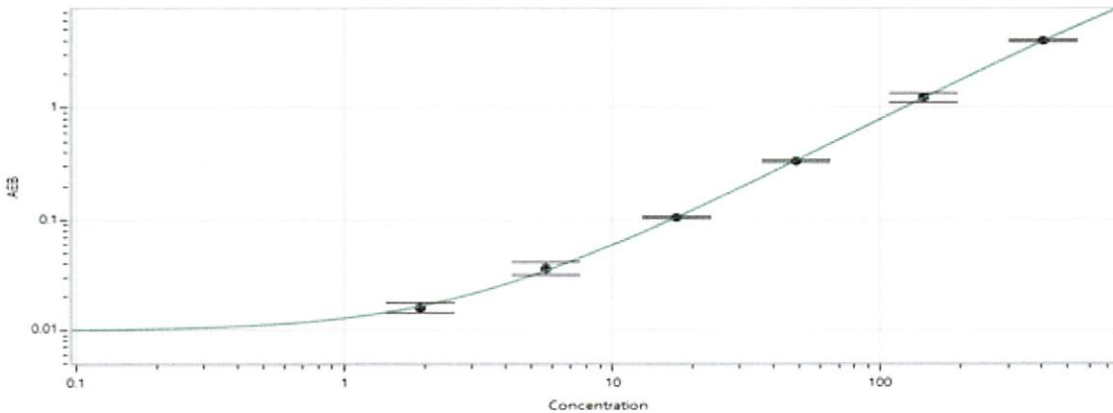
¹ Reagents are not Kit Lot Specific

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

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	333804	41.8	35.3 - 52.9
Control 2	333804	1135	869 - 1303
Panel 1	301702	14.8	12.7 - 20.9
Panel 2	301702	14.0	12.3 - 20.9
Panel 3	309704	416	331 - 645
Panel 4	301702	99.9	57.9 - 125

* 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	1.91
C	5.69
D	17.3
E	48.2
F	144
G	406

Review/Approval

Cody Carlson

QC Manager

Name

Title

Cody Carlson 16 Jul 2024
Signature/Date