

Product Number:	104111
Lot Number:	504132
Expiration Date:	21-Sep-25
Platform(s):	SR-X and HD-1/HD-X



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

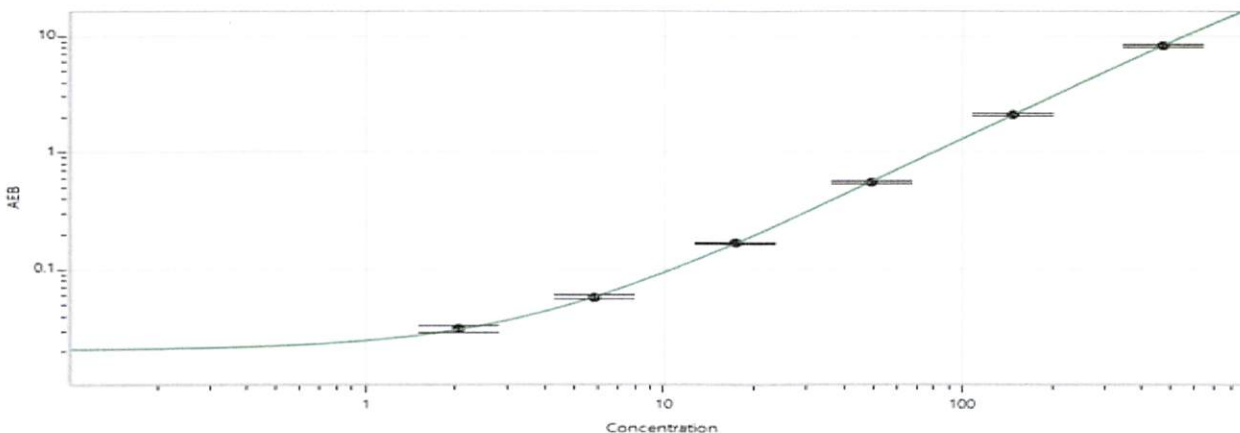
Component	Part Number	Lot Number
Bead Reagent	104098	408109
Detector Reagent	104099	408607
SBG Reagent	104100	408209
Sample Diluent	104101	408112
Calibrators	104106	409203
Control 1	104107	409204
Control 2	104108	409204
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	409204	39.6	32.2 - 48.3
Control 2	409204	1054	840 - 1260
Panel 1	301702	14.2	12.7 - 20.9
Panel 2	301702	14.1	12.3 - 20.9
Panel 3	309704	406	331 - 645
Panel 4	406103	194	157 - 236

* 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	2.04
C	5.85
D	17.3
E	49.5
F	148
G	470

Review/Approval

Cody Carlson

QC Manager

Name

Title

Cody Carlson 25 Jun 2024
Signature/Date