

<b>Product Number:</b>	100860
<b>Lot Number:</b>	504176
<b>Expiration:</b>	25-Jan-2025
<b>Platform(s):</b>	HD-1/HD-X



Component	Part Number	Lot Number
Bead Reagent	100865	320710
Detector Reagent	100862	321208
SBG Reagent	100868	323008
Sample Diluent	100870	320605
Calibrators	101111	320902
Control 1	100875	320903
Control 2	100876	320903
RGP Reagent	103159	N/A <sup>1</sup>

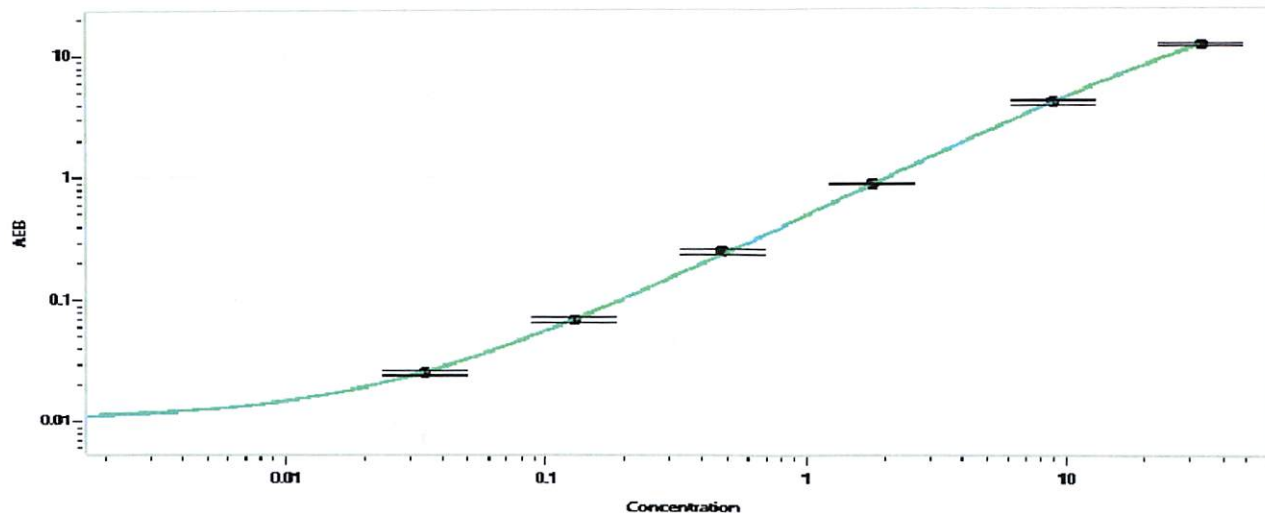
<sup>1</sup> Reagents are not Kit Lot Specific

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

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	320903	6.04	5.23 - 8.10
Control 2	320903	0.097	0.075 - 0.127
Panel 1	306704	0.794	0.772 - 1.16
Panel 2	306704	3.94	2.98 - 7.27
Panel 3	306704	15.7	10.1 - 25.4

\* 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	0.034
C	0.128
D	0.481
E	1.79
F	8.98
G	32.9

Review/Approval

Cody Carlson

QC Manager

*[Signature]* 06 MAR 2024

Name

Title

Signature/Date