

Product Number:	101599
Lot Number:	502651
Expiration:	7-Dec-2021
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	101601	034209
Detector Reagent	102012	103003
SBG Reagent	101602	100902
Sample Diluent	101603	034602
Calibrator Diluent	102014	031416
Calibrator Concentrate	100778	001414
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific



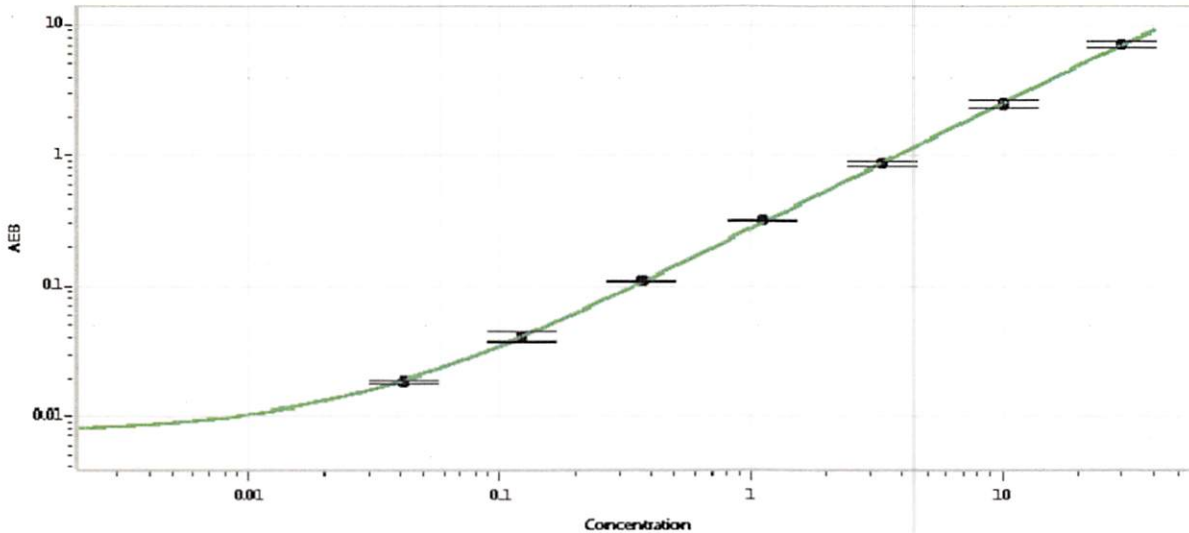
Data below represents results generated on the Simoa™ HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-012518-02	0.388	0.352-0.537
Control 2	MF-012518-02	21.9	14.5-28.3
Panel 3	034908	9.08	7.63-12.2
Panel 4	034908	35.2	24.7-37.0

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

Calibrator Concentrate Value Assignment (pg/mL)	706
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Junpil Park

QC Supervisor

Name

Title

Signature/Date

[Handwritten Signature] 25 Feb 2021