

Product Number:	102860
Lot Number:	502925
Expiration:	29-Sep-2021
Platform(s):	HD-1/HD-X

Component	Part Number	Lot Number
Bead Reagent	102852	113202
Detector Reagent	102854	108802
SBG Reagent	102855	108803
Sample Diluent	102856	108804
Calibrator Diluent	102859	108805
Calibrator Concentrate	102857	002401
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific



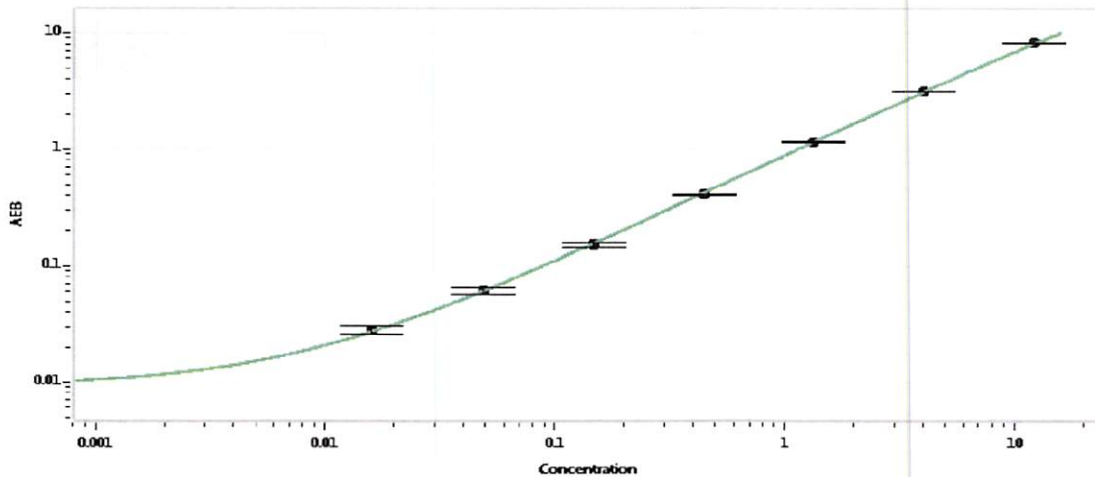
Data below represents results generated on the Simoa™ HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	002401	0.229	0.180-0.300
Control 2	002401	4.09	3.60-6.00
Panel 1	033802	0.148	0.118-0.249
Panel 2	033802	1.95	1.64-2.65
Panel 3	033802	9.12	7.98-12.0

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

Calibrator Concentrate Value Assignment (pg/mL)	240
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.016
C	0.049
D	0.148
E	0.444
F	1.33
G	4.00
H	12.0

Review/Approval

Junpil Park

QC Supervisor

 04 Jun 2021

Name

Title

Signature/Date