

Product Number:	101599
Lot Number:	502794
Expiration:	4-Mar-2022
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	101601	106404
Detector Reagent	102012	106605
SBG Reagent	101602	106601
Sample Diluent	103886	106602
Calibrator Diluent	102014	106316
Calibrator Concentrate	100778	106707
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific



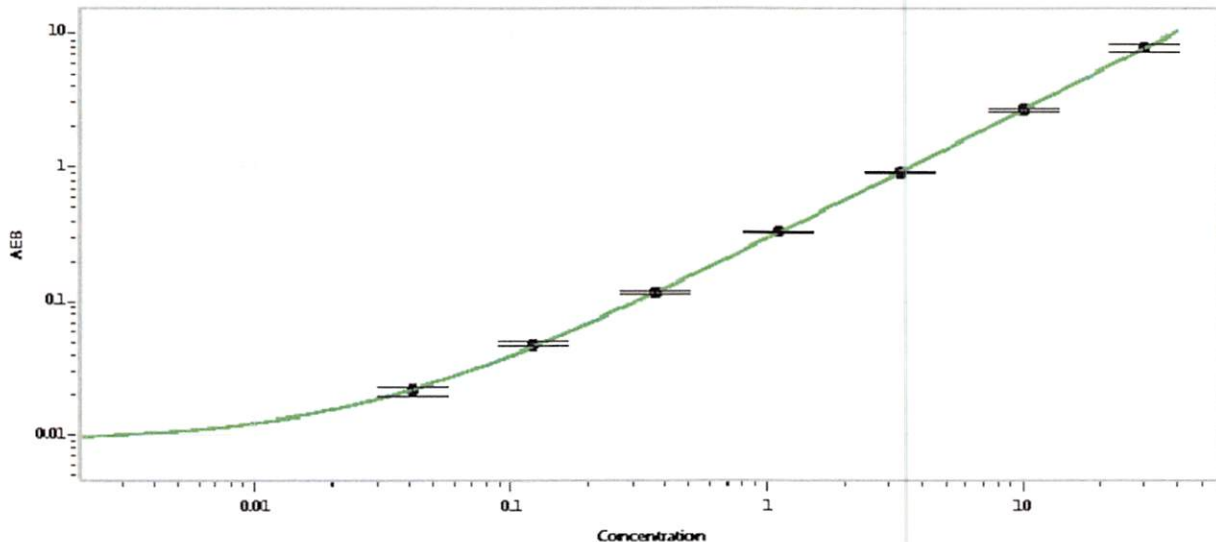
Data below represents results generated on the Simoa™ HD-X Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	108214	0.259	0.223-0.374
Control 2	109503	42.2	29.4-46.0
Panel 3	034908	11.3	7.76-12.7
Panel 4	034908	40.7	20.9-45.7

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

Calibrator Concentrate Value Assignment (pg/mL)	1004
---	------

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

[Signature] 15 Apr 2021

Name

Title

Signature/Date