

Product Number:	100794
Lot Number:	502860
Expiration:	19-Nov-2021
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Reagent	100781	109818
Detector Reagent	100782	110320
SBG Reagent	100783	110003
Sample Diluent	103202	107722
Calibrator Diluent	100793	108127
Calibrator Concentrate	100780	932311
SBG Diluent (SRX)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific



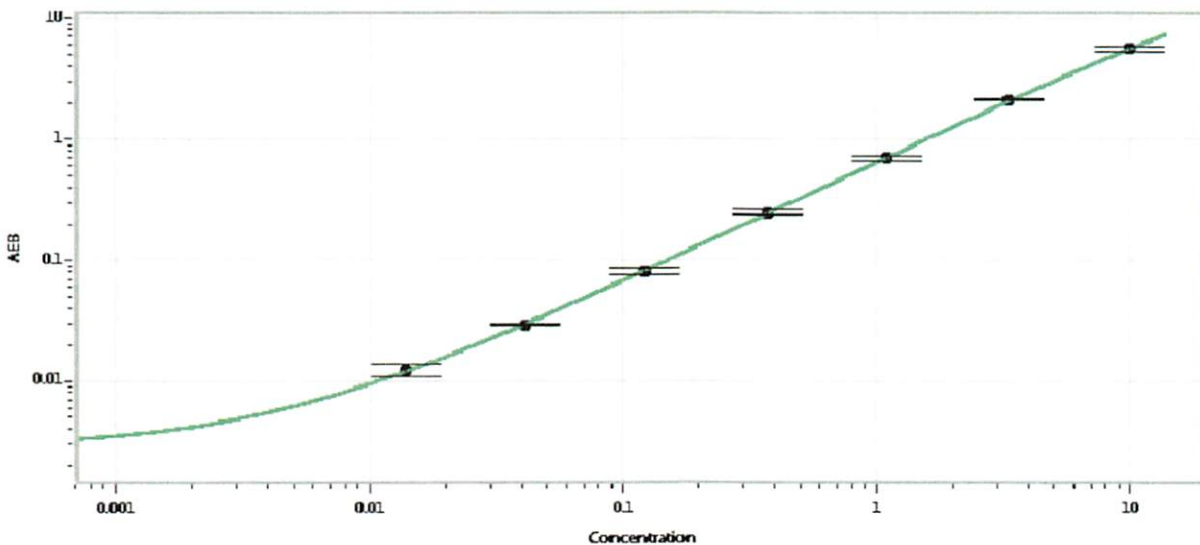
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	920002	0.229	0.215-0.356
Control 2	920002	3.09	2.88-4.31
Panel 1	100801	2.52	2.22-3.45
Panel 2	100801	16.8	14.5-21.8
Panel 3	100801	1.80	1.15-2.98

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

Calibrator Concentrate Value Assignment (pg/mL)	134
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.014
C	0.041
D	0.123
E	0.370
F	1.11
G	3.33
H	10.0

Review/Approval

Junpil Park

QC Supervisor

[Handwritten Signature] 26 Apr 2021

Name

Title

Signature/Date