

Product Number:	101664
Lot Number:	502914
Expiration:	2-Apr-2022
Platform(s):	HD-1/HD-X

Component	Part Number	Lot Number
Bead Reagent	101666	112405
Detector Reagent	101665	112407
SBG Reagent	101667	112408
Sample Diluent	101668	109218
Calibrator Diluent	101669	112409
Calibrator Concentrate	103217	106408
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific



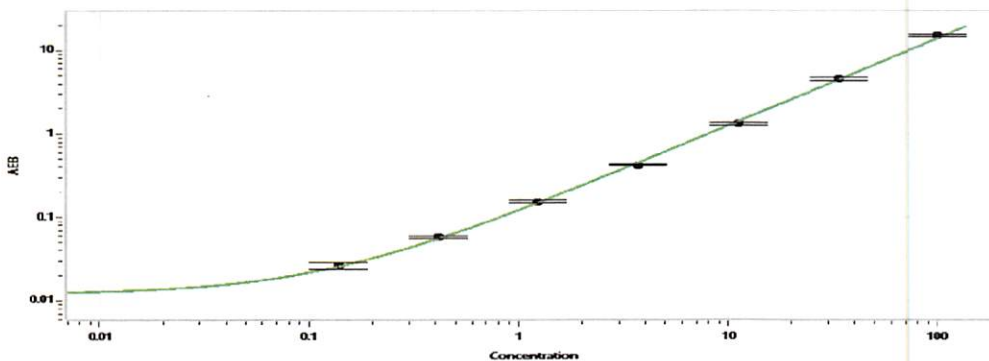
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	001402	0.674	0.284-1.19
Control 2	001402	32.2	17.8-59.8
Panel 1	934301	5.55	3.30-9.85
Panel 2	934302	21.2	12.3-40.5

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

Calibrator Concentrate Value Assignment (pg/mL)	1674
---	------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.137
C	0.412
D	1.23
E	3.70
F	11.1
G	33.3
H	100

Review/Approval

Junpil Park
Name

QC Supervisor
Title

[Signature] 20 May 2021
Signature/Date