

Product Number:	101664
Lot Number:	502928
Expiration:	10-May-2022
Platform(s):	HD-1/HD-X

Component	Part Number	Lot Number
Bead Reagent	101666	113009
Detector Reagent	101665	113111
SBG Reagent	101667	113112
Sample Diluent	101668	113113
Calibrator Diluent	101669	113114
Calibrator Concentrate	103217	113306
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific



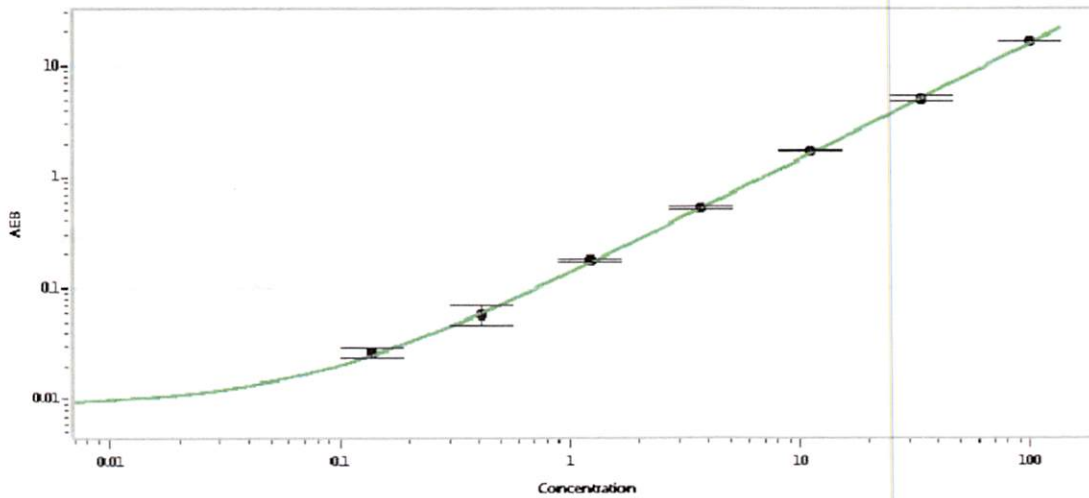
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	001402	0.689	0.316-1.14
Control 2	001402	33.5	18.0-57.8
Panel 1	934301	6.45	3.30-9.85
Panel 2	934302	24.5	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)	1865
---	------

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

QC Supervisor

[Signature] 16 Jan 2021

Name

Title

Signature/Date