

Product Number:	101664
Lot Number:	502904
Expiration:	23-Sep-2021
Platform(s):	HD-1/HD-X



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

Component	Part Number	Lot Number
Bead Reagent	101666	105513
Detector Reagent	101665	111502
SBG Reagent	101667	105609
Sample Diluent	101668	105611
Calibrator Diluent	101669	026723
Calibrator Concentrate	103217	106408
RGP Reagent	103159	N/A ¹

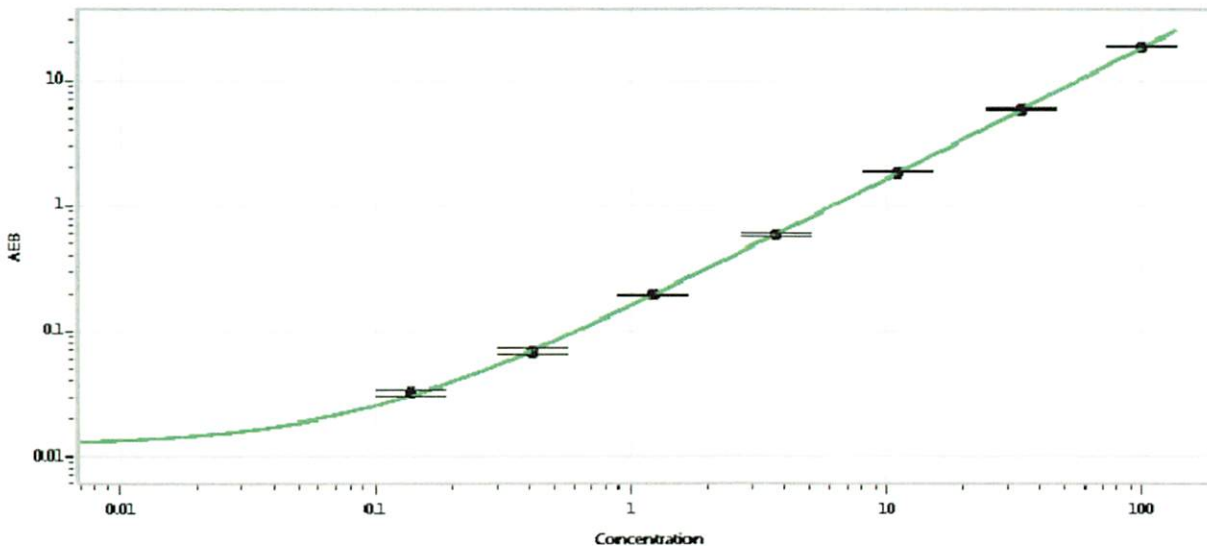
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	001402	0.622	0.261-1.30
Control 2	001402	33.1	19.2-62.8
Panel 1	934301	6.28	3.41-10.1
Panel 2	934302	22.8	13.2-41.3

* 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/Date
03 May 2021