

Product Number:	101622
Lot Number:	502873
Expiration:	3-Feb-2022
Platform(s):	HD-1 / HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101624	110513
Detector Reagent	101623	105102
SBG Reagent	101625	106418
Sample Diluent	103732	103416
Calibrator Diluent	101627	105012
Calibrator Concentrate	100652	105604
RGP Reagent	103159	N/A <sup>1</sup>

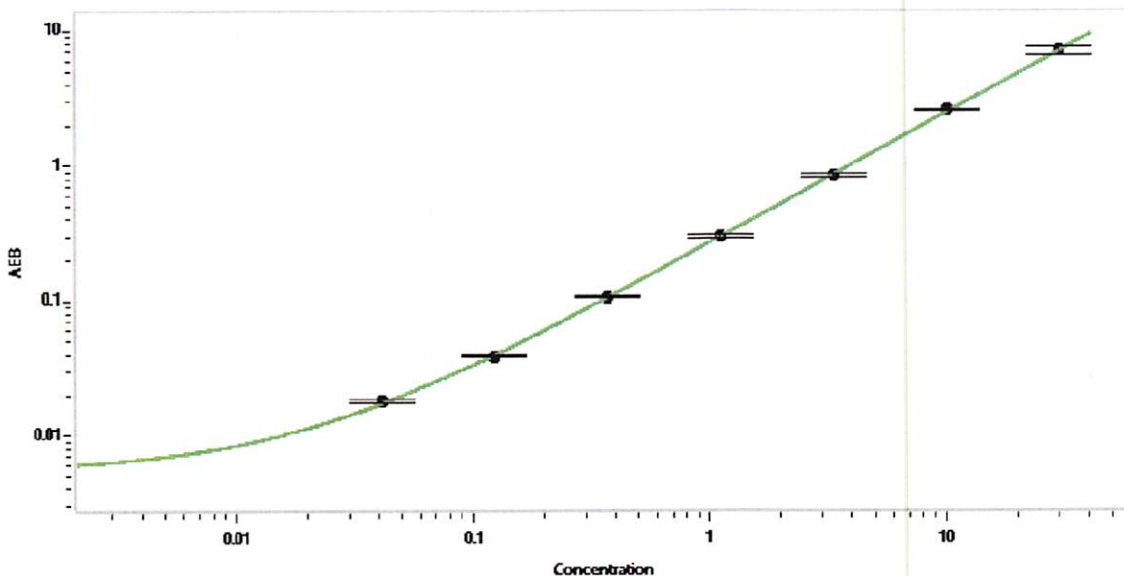
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	108406	1.20	1.12-2.04
Control 2	108406	23.4	17.2-25.9
Panel 1	029301	45.6	39.8-59.7
Panel 2	029401	4.68	4.35-6.53
Panel 3	029301	1.28	1.22-1.82

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

Calibrator Concentrate Value Assignment (pg/mL)	601
---	-----

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  
Name

QC Supervisor  
Title

 17 May 2021  
Signature/Date