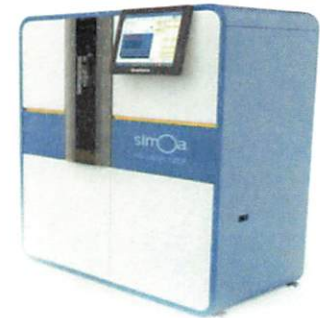


Product Number:	101622
Lot Number:	503042
Expiration:	20-Jun-2022
Platform(s):	HD-1 / HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101624	119405
Detector Reagent	101623	119412
SBG Reagent	101625	119413
Sample Diluent	103732	117118
Calibrator Diluent	101627	119414
Calibrator Concentrate	100652	119411
RGP Reagent	103159	N/A <sup>1</sup>

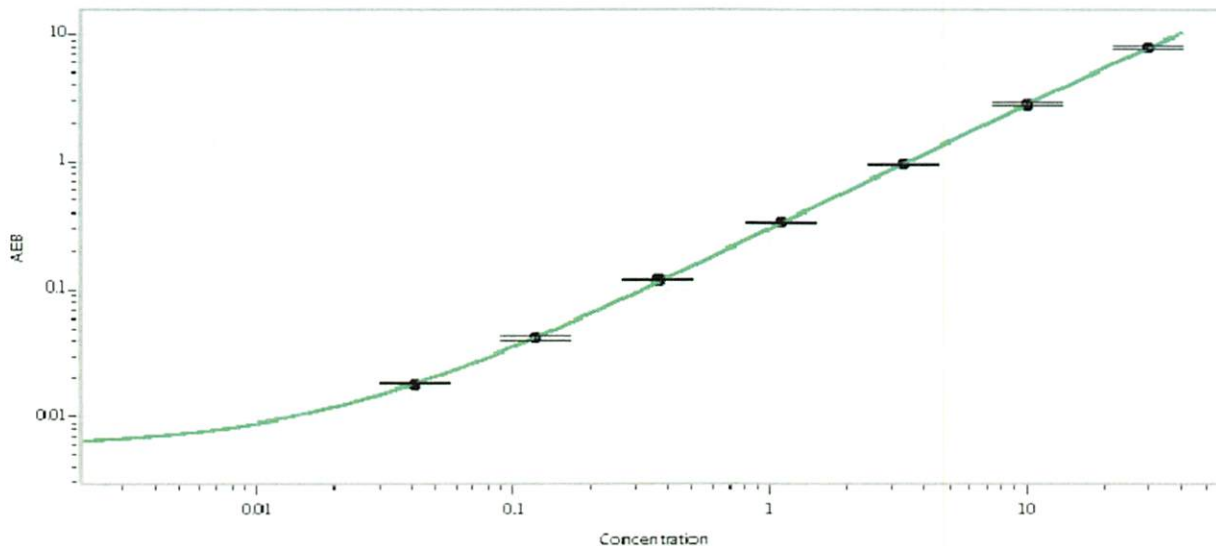
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	108406	1.63	1.19-1.99
Control 2	108406	26.1	16.4-28.8
Panel 1	029301	57.5	39.8-59.7
Panel 2	029401	6.36	4.35-6.53
Panel 3	029301	1.74	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)	763
---	-----

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

QC Supervisor

*[Signature]* 22 Jul 2021

Name

Title

Signature/Date