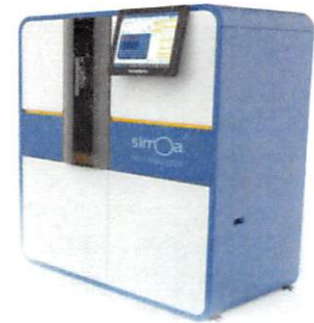


Product Number:	101599
Lot Number:	503033
Expiration:	29-Jun-2022
Platform(s):	HD-1/HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101601	118901
Detector Reagent	102012	119001
SBG Reagent	101602	119002
Sample Diluent	103886	119003
Calibrator Diluent	102014	118022
Calibrator Concentrate	100778	119004
RGP Reagent	103159	N/A ¹

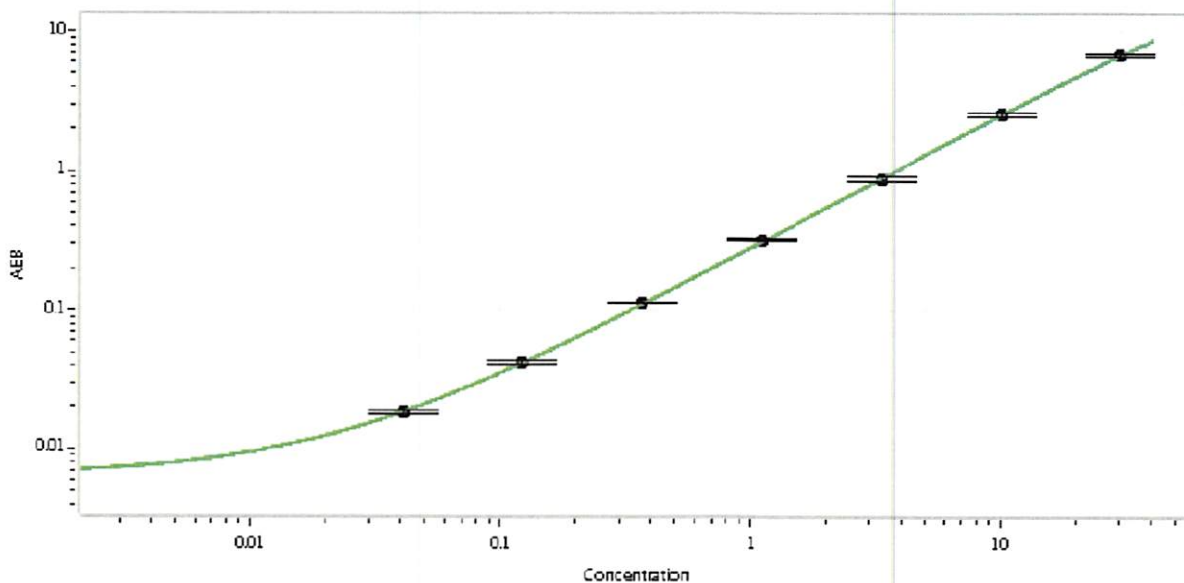
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	108214	0.194	0.146-0.396
Control 2	109503	31.6	28.7-47.3
Panel 3	034908	9.07	8.12-12.5
Panel 4	034908	32.5	22.4-47.5

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

Calibrator Concentrate Value Assignment (pg/mL)	555
---	-----

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

23 Jul 2021
Signature/Date