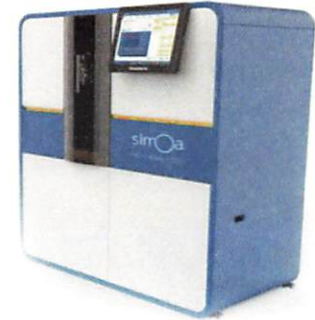


Product Number:	102329
Lot Number:	503145
Expiration:	1-Jun-2022
Platform(s):	HD-1/HD-X



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

Component	Part Number	Lot Number
Bead Reagent	102322	126308
Detector Reagent	102324	126309
SBG Reagent	102325	126311
Sample Diluent	102326	115233
Calibrator Diluent	102327	121312
Calibrator Concentrate	100849	126310
RGP Reagent	103159	N/A ¹

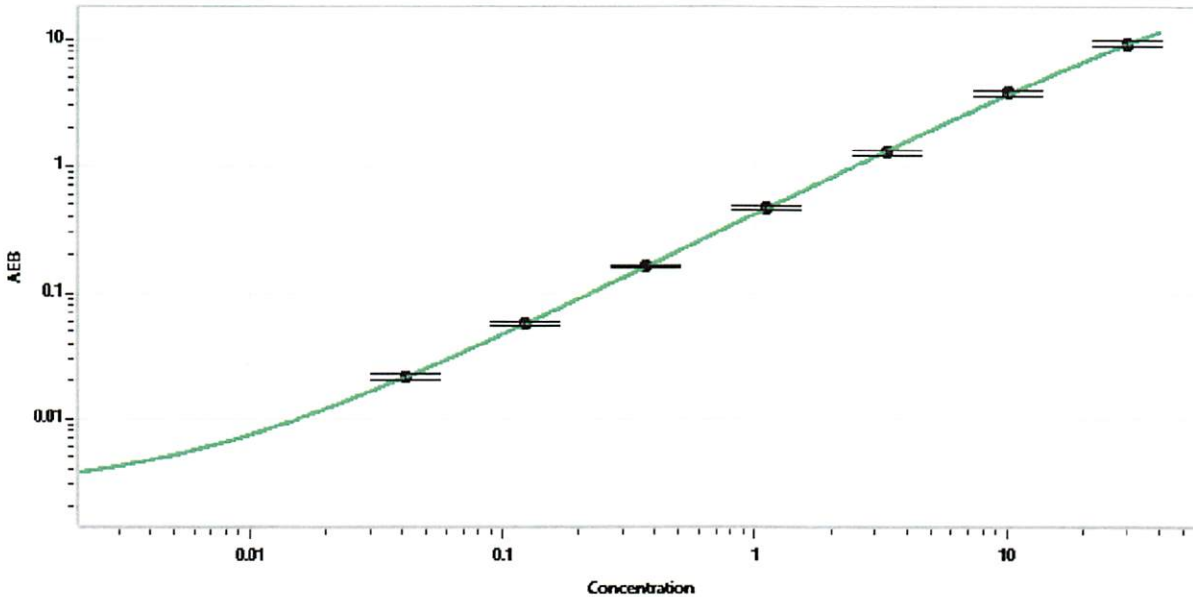
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	023702	0.201	0.114-0.249
Control 2	023702	23.0	17.6-33.7
Panel 1	027607	0.145	0.050-0.176
Panel 2	027607	0.154	0.071-0.213
Panel 3	027607	11.0	7.45-18.2

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

Calibrator Concentrate Value Assignment (pg/mL)	519
-------------------------------------------------	-----

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] 14Oct-2021

Name

Title

Signature/Date