

Product Number:	102329
Lot Number:	502903
Expiration:	15-Feb-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	104715
Detector Reagent	102324	104803
SBG Reagent	102325	111016
Sample Diluent	102326	104618
Calibrator Diluent	102327	104723
Calibrator Concentrate	100849	108120
RGP Reagent	103159	N/A <sup>1</sup>

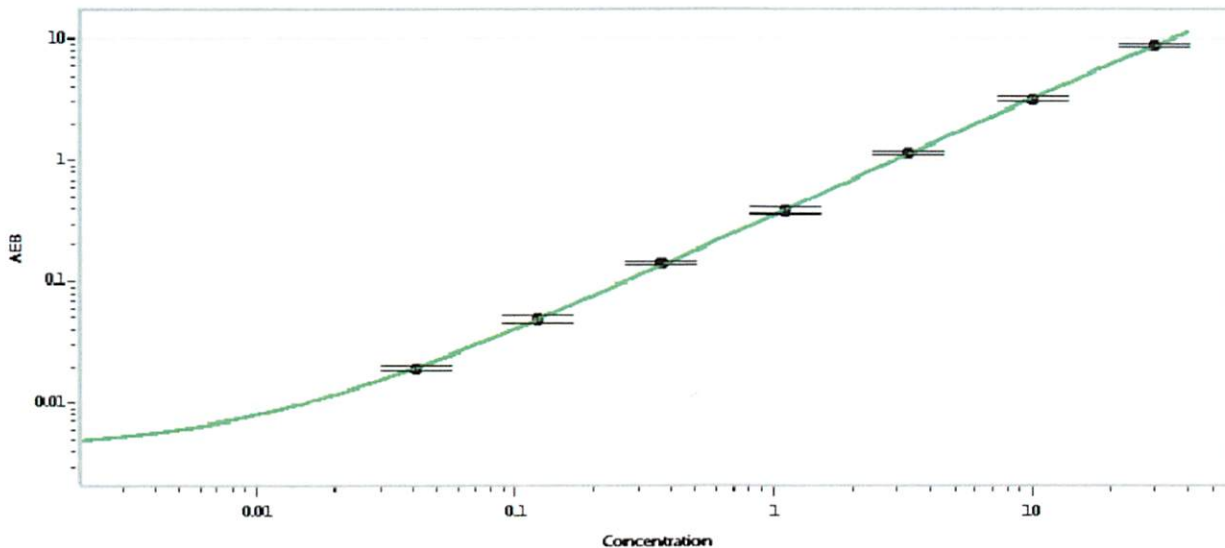
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	023702	0.192	0.119-0.229
Control 2	023702	26.9	18.8-30.7
Panel 1	027607	0.108	0.037-0.142
Panel 2	027607	0.106	0.063-0.168
Panel 3	027607	14.4	7.66-17.0

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

Calibrator Concentrate Value Assignment (pg/mL)	5585
---	------

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

 21 May 2021  
Signature/Date