

Product Number:	101599
Lot Number:	503417
Expiration:	16-Feb-2023
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	101601	213305
Detector Reagent	102012	213309
SBG Reagent	101602	213310
Sample Diluent	103886	213704
Calibrator Diluent	102014	210116
Calibrator Concentrate	100778	204206
RGP Reagent	103159	N/A <sup>1</sup>

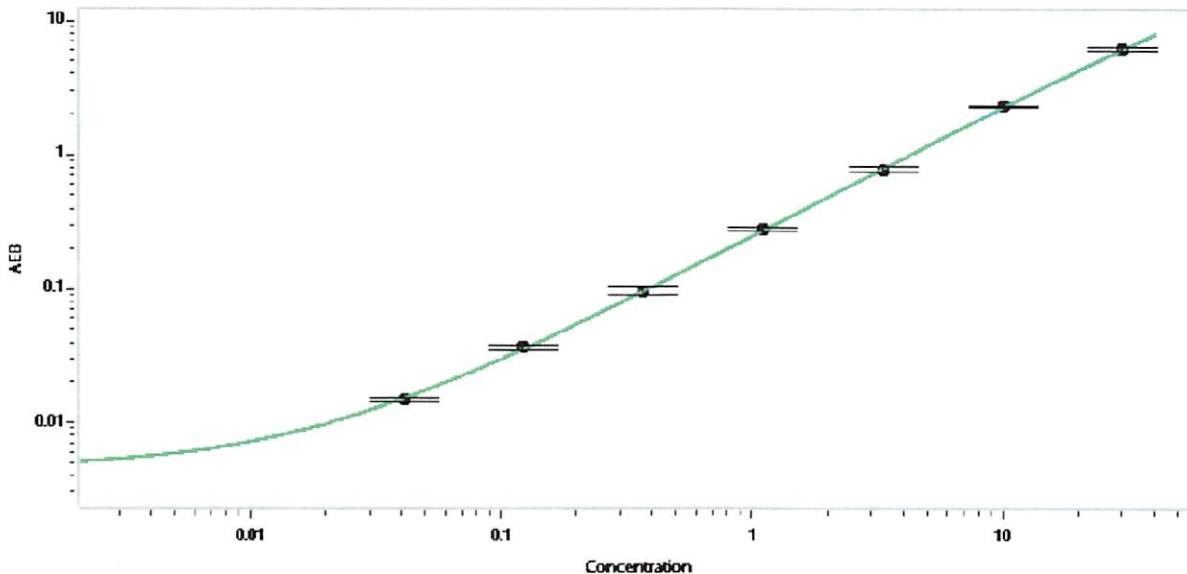
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	204206	23.1	18.0-30.0
Control 2	204206	0.471	0.375-0.625
Panel 3	034908	10.7	8.12-12.5
Panel 4	034908	34.7	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)	574
---	-----

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

  
 Name
 

 Director, QC  
 Title
 

  
 Signature/Date
 

 16 Jun 2022