

Product Number:	101599
Lot Number:	502997
Expiration:	27-Apr-2022
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	113805
Detector Reagent	102012	114807
SBG Reagent	101602	113816
Sample Diluent	103886	116513
Calibrator Diluent	102014	111726
Calibrator Concentrate	100778	113904
RGP Reagent	103159	N/A <sup>1</sup>

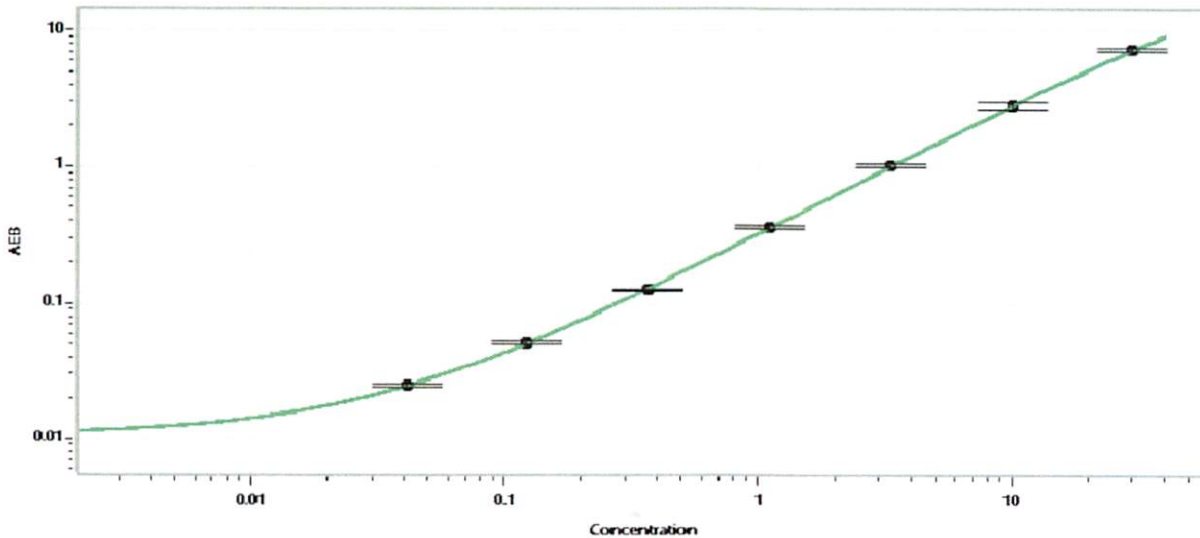
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	108214	0.219	0.146-0.396
Control 2	109503	36.6	28.7-47.3
Panel 3	034908	8.47	7.91-12.5
Panel 4	034908	30.5	22.3-45.6

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

Calibrator Concentrate Value Assignment (pg/mL)	559
---	-----

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

  
Signature/Date

24 Jun 2021