

Product Number:	102860
Lot Number:	503328
Expiration:	6-Aug-2022
Platform(s):	HD-1/HD-X

Component	Part Number	Lot Number
Bead Reagent	102852	203701
Detector Reagent	102854	203813
SBG Reagent	102855	203814
Sample Diluent	102856	203817
Calibrator Diluent	102859	203816
Calibrator Concentrate	102857	203913
RGP Reagent	103159	N/A <sup>1</sup>



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

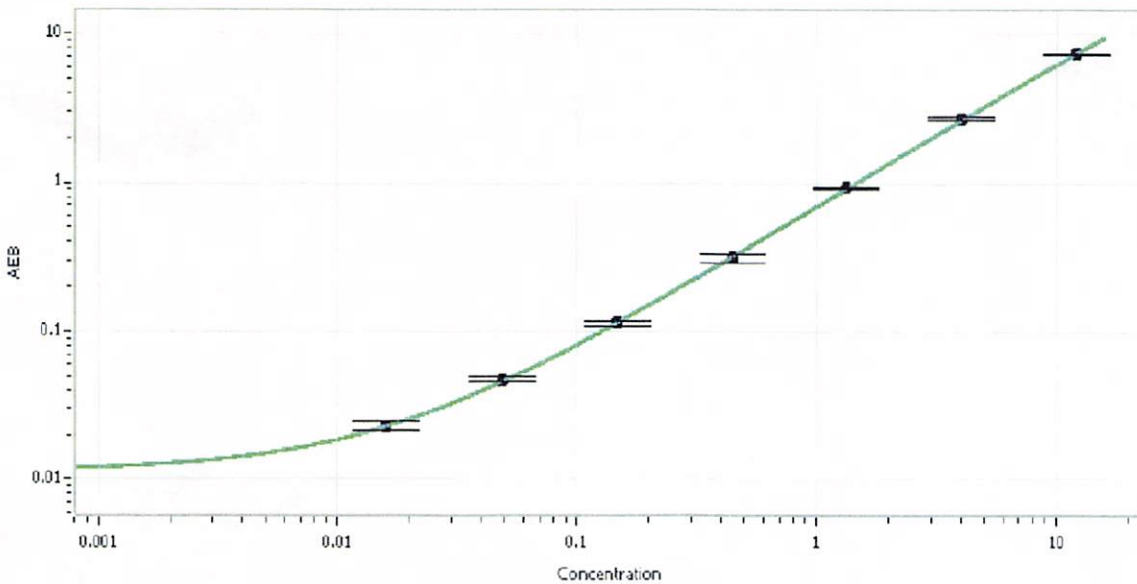
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	121506	0.258	0.229 - 0.344
Control 2	121506	4.83	3.93 - 5.90
Panel 1	033802	0.132	0.109 - 0.242
Panel 2	033802	1.97	1.70 - 2.61
Panel 3	033802	9.45	8.09 - 12.1

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

Calibrator Concentrate Value Assignment (pg/mL)	307
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.016
C	0.049
D	0.148
E	0.444
F	1.33
G	4.00
H	12.0

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel 24 Feb 2022  
Signature/Date