

Product Number:	101622
Lot Number:	503389
Expiration:	17-Sep-2022
Platform(s):	HD-1 / HD-X & SR-X



Component	Part Number	Lot Number
Bead Reagent	101624	126608
Detector Reagent	101623	203204
SBG Reagent	101625	127007
Sample Diluent	103732	126007
Calibrator Diluent	101627	204614
Calibrator Concentrate	100652	203712
RGP Reagent	103159	N/A <sup>1</sup>

Data below represents results generated on the Simoa® HD-1 / HD-X & SR-X Analyzer.

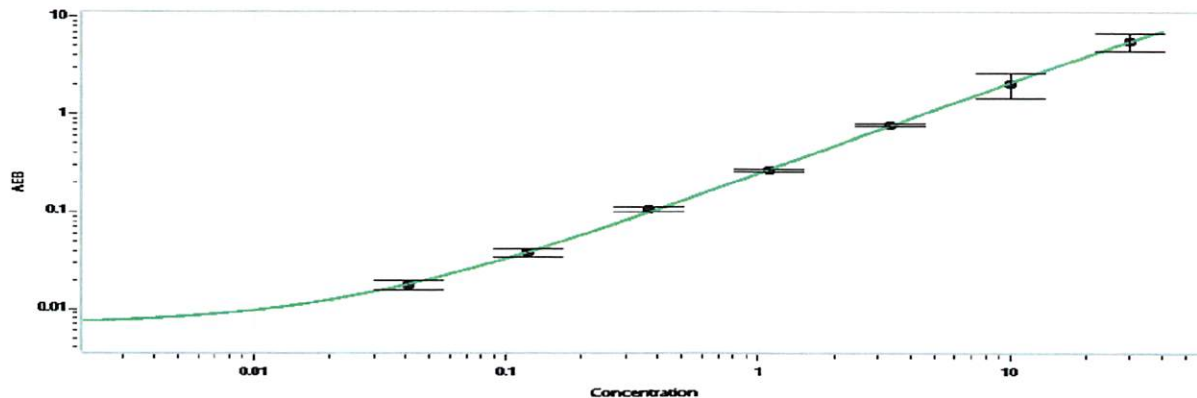
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	203712	19.4	15.0-25.0
Control 2	203712	2.15	1.50-2.50
Panel 1	029301	46.0	39.8-59.7
Panel 2	029401	4.78	4.35-6.53
Panel 3	029301	1.23	1.22-1.82

\* 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

*Shivani Goel*  
Shivani Goel

QC Manager

*Shivani Goel 12 May 2022*  
Signature/Date