

Product Number:	101622
Lot Number:	503299
Expiration:	5-Aug-2022
Platform(s):	HD-1 / HD-X & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101624	126608
Detector Reagent	101623	203204
SBG Reagent	101625	127007
Sample Diluent	103732	126007
Calibrator Diluent	101627	121720
Calibrator Concentrate	100652	203712
RGP Reagent	103159	N/A <sup>1</sup>

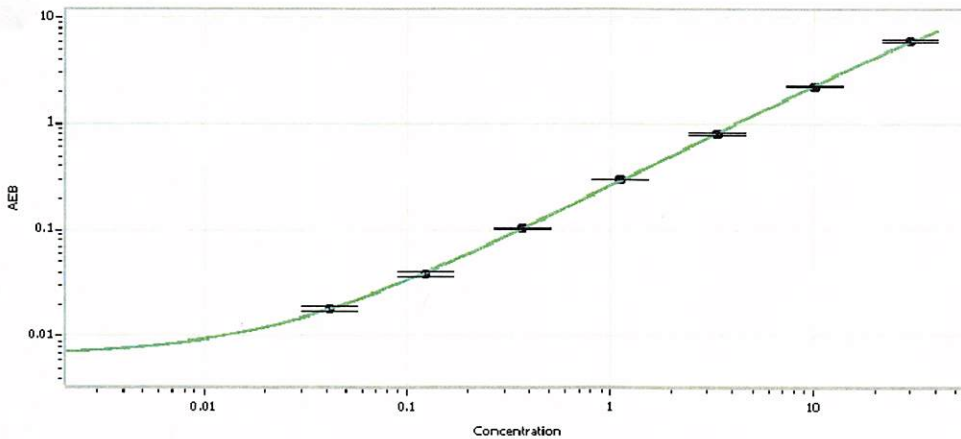
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	108406	1.49	1.23-1.92
Control 2	108406	25.0	17.5-28.9
Panel 1	029301	52.2	39.8-59.7
Panel 2	029401	5.71	4.35-6.53
Panel 3	029301	1.53	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*  
Name

QC  
Title

*Shivani Goel 24Feb2022*  
Signature/Date