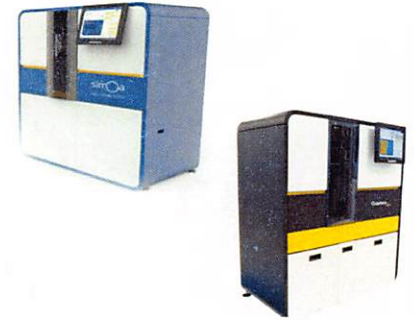


Product Number:	101672
Lot Number:	503362
Expiration:	7-Oct-2022
Platform(s):	HD-1/HD-X



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

Component	Part Number	Lot Number
Bead Reagent	101674	129704
Detector Reagent	101673	129803
SBG Reagent	101675	129907
Sample Diluent	101676	133927
Calibrator Diluent	101677	128016
Calibrator Concentrate	101734	130203
RGP Reagent	103159	N/A <sup>1</sup>

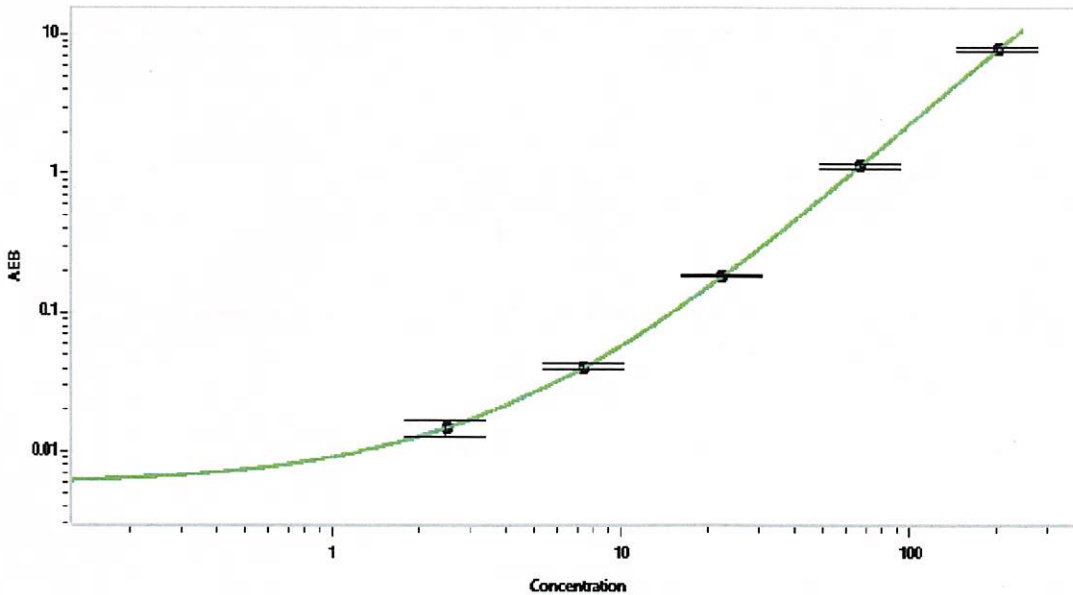
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	126904	19.6	13.2-23.9
Control 2	126904	399	308-484
Panel 1	024003	120	94-149
Panel 2	024003	248	198-326

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

Calibrator Concentrate Value Assignment (pg/mL)	4295
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.47
C	7.41
D	22.2
E	66.7
F	200

Review/Approval

Matthew Smart  
Name

QC Director  
Title

*MSA* 24 Mar 2022  
Signature/Date