

Product Number:	103277
Lot Number:	502229
Expiration:	21-Nov-2020
Platform(s):	HD-1/ HDX



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

Component	Part Number	Lot Number
Bead Reagent	103265	001706
Detector Reagent	103268	003509
SBG Reagent	103269	000713
Sample Diluent	103270	000805
Calibrator Diluent	103266	000806
Calibrator Concentrate	103272	832501
RGP Reagent	103159	N/A <sup>1</sup>

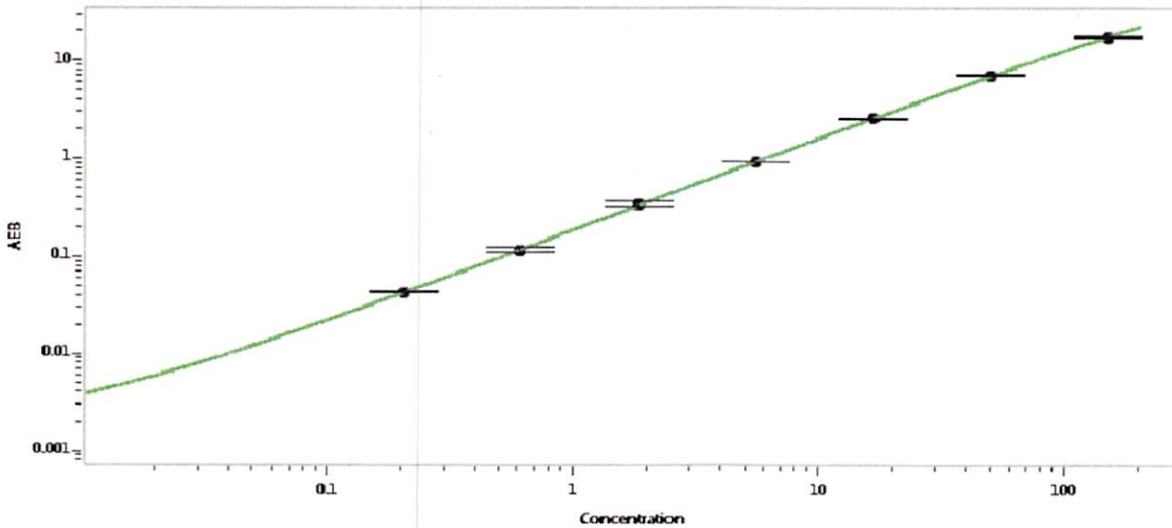
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	824302	4.76	4.16-6.24
Control 2	824302	137	111-167
Panel 1	823501	2.11	1.77-3.02
Panel 2	823502	28.9	21.7-37.1
Panel 3	823503	98.0	72.5-109

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

Calibrator Concentrate Value Assignment (pg/mL)	5550
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.206
C	0.617
D	1.85
E	5.56
F	16.7
G	50.0
H	150

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P Stadler</i>	07 Feb 2020
	Name	Title	Signature/Date	