

Product Number:	103277
Lot Number:	501523
Expiration:	20-Nov-2019
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	103265	832404
Detector Reagent	103268	832505
SBG Reagent	103269	832504
Sample Diluent	103270	832402
Calibrator Diluent	103266	832403
Calibrator Concentrate	103272	832501
RGP Reagent	103159	N/A ¹

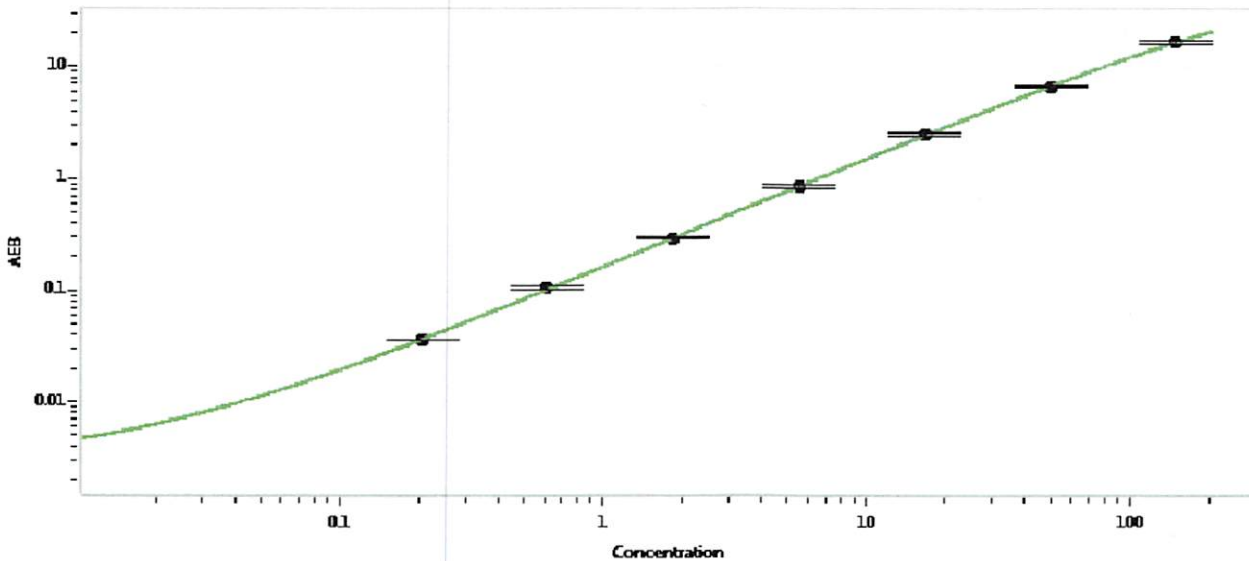
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	824302	5.01	4.22-6.33
Control 2	824302	131	111-167
Panel 1	823501	2.44	1.85-3.04
Panel 2	823502	30.4	21.3-37.8
Panel 3	823503	92.9	71.6-107

* 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	Associate Director, QC	<i>P. Stadler</i>	11/6/2019
	Name	Title	Signature/Date	