

Product Number:	101622
Lot Number:	501694
Expiration:	29-Jan-2020
Platform(s):	HD-1 & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101624	905811
Detector Reagent	101623	905806
SBG Reagent	101625	905807
Sample Diluent	103203	902930
Calibrator Diluent	101627	904416
Calibrator Concentrate	100652	906001
RGP Reagent	103159	N/A ¹

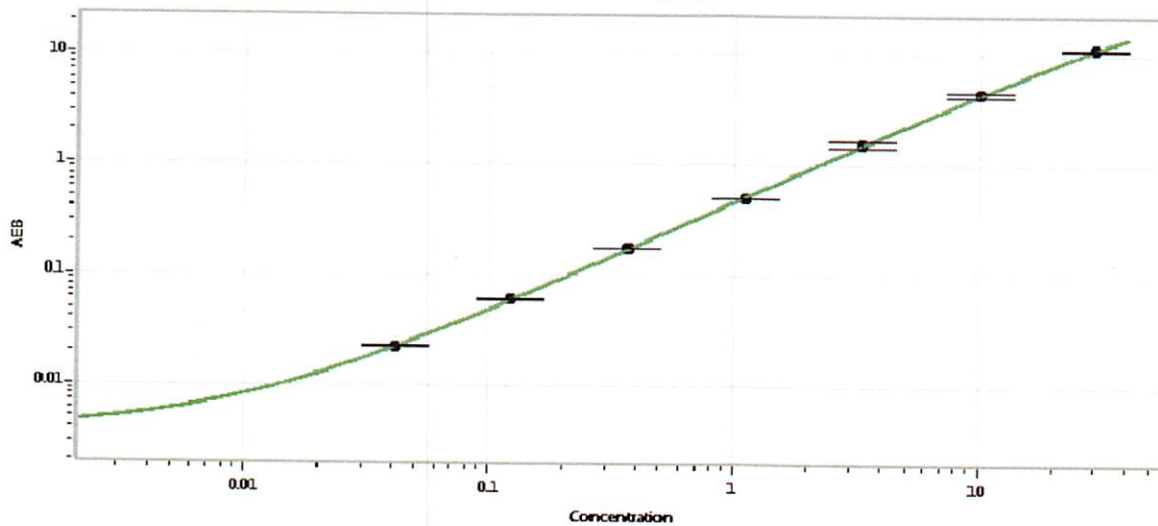
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	901702	1.23	0.905-1.36
Control 2	901702	19.6	14.5-21.7
Panel 1	MF-050918-01	55.3	41.3-66.4
Panel 2	MF-050918-02	7.50	5.08-8.55
Panel 3	MF-050918-03	1.04	0.635-1.22

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

Calibrator Concentrate Value Assignment (pg/mL)	864
---	-----

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	Patrizia Stadler	Associate Director, QC	<i>P. Stadler</i>	07Mar2019
	Name	Title	Signature/Date	