

Product Number:	101664
Lot Number:	501936
Expiration:	29-Jan-2020
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	101666	903607
Detector Reagent	101665	903608
SBG Reagent	101667	903104
Sample Diluent	101668	902910
Calibrator Diluent	101669	902911
Calibrator Concentrate	103217	834607
RGP Reagent	103159	N/A ¹

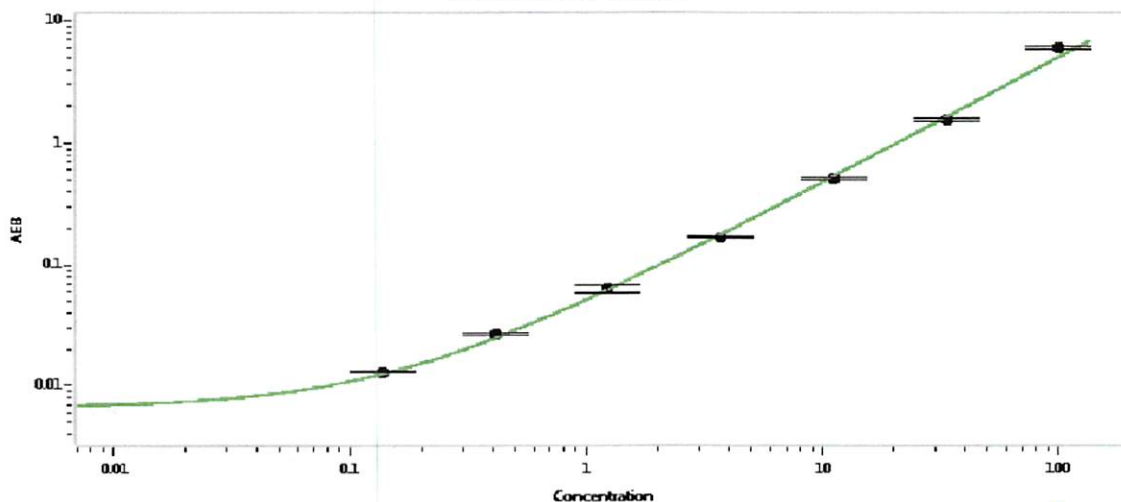
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	52520	1.26	0.740-1.33
Control 2	52520	35.3	27.1-40.6
Panel 1	MF-082015-01	1.39	0.942-1.53
Panel 2	MF-081415-02	10.2	7.46-11.2
Panel 3	MF-081415-03	12.1	9.20-15.0
Panel 4	MF-090315-09	45.9	27.6-66.8

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

Calibrator Concentrate Value Assignment (pg/mL)	1427
---	------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.137
C	0.412
D	1.23
E	3.70
F	11.1
G	33.3
H	100

Review/Approval

Daniel Ritchie QC Senior Analyst
Name Title

[Signature] 1/23/2019
Signature/Date