

Product Number:	101580
Lot Number:	501756
Expiration:	14-Mar-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	101582	908804
Detector Reagent	101581	908511
SBG Reagent	101583	908407
Sample Diluent	103201	907309
Calibrator Diluent	101585	909205
Calibrator Concentrate	100655	834101
RGP Reagent	103159	N/A <sup>1</sup>

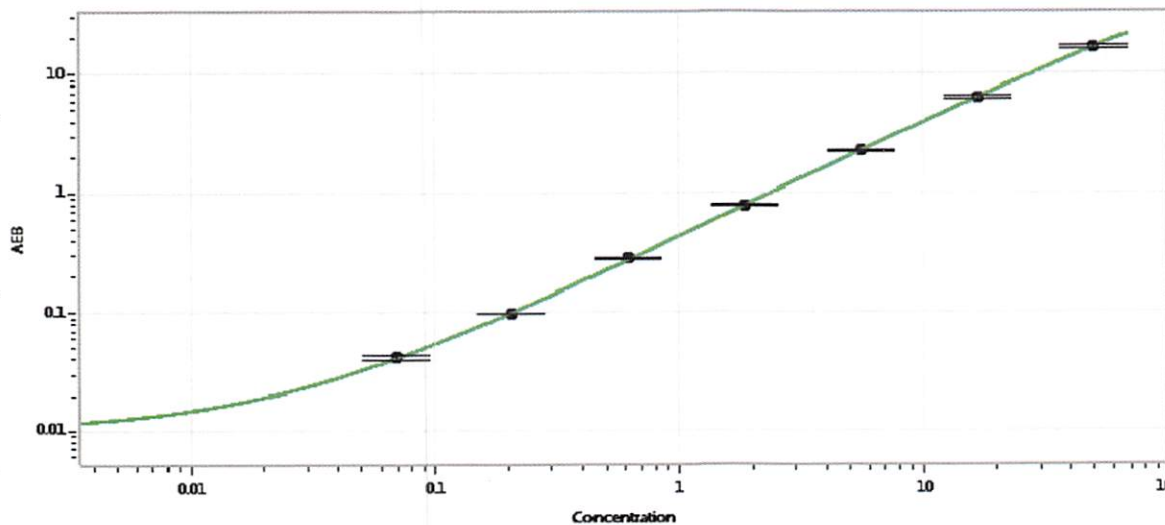
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-042817-01	2.77	2.05-3.08
Control 2	MF-051617-02	50.2	35.8-53.7
Panel 1	MF-051017-03	28.0	20.9-32.4
Panel 2	MF-032217-11	1.56	1.33-2.15
Panel 3	MF-033018-06	2.59	1.68-3.50

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

Calibrator Concentrate Value Assignment (pg/mL)	1155 pg/mL
---	------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.069
C	0.206
D	0.617
E	1.85
F	5.56
G	16.7
H	50.0

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i> 05 Apr 2019
	Name	Title	