

Product Number:	101672
Lot Number:	501177
Expiration:	19-Sep-2018
Platform(s):	HD-1



Component	Part Number	Lot Number
Bead Reagent	101674	80803
Detector Reagent	101673	80782
SBG Reagent	101675	80805
Sample Diluent	101676	73499
Calibrator Diluent	101677	80474
Calibrator Concentrate	101734	73205
RGP Reagent	103159	N/A <sup>1</sup>

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

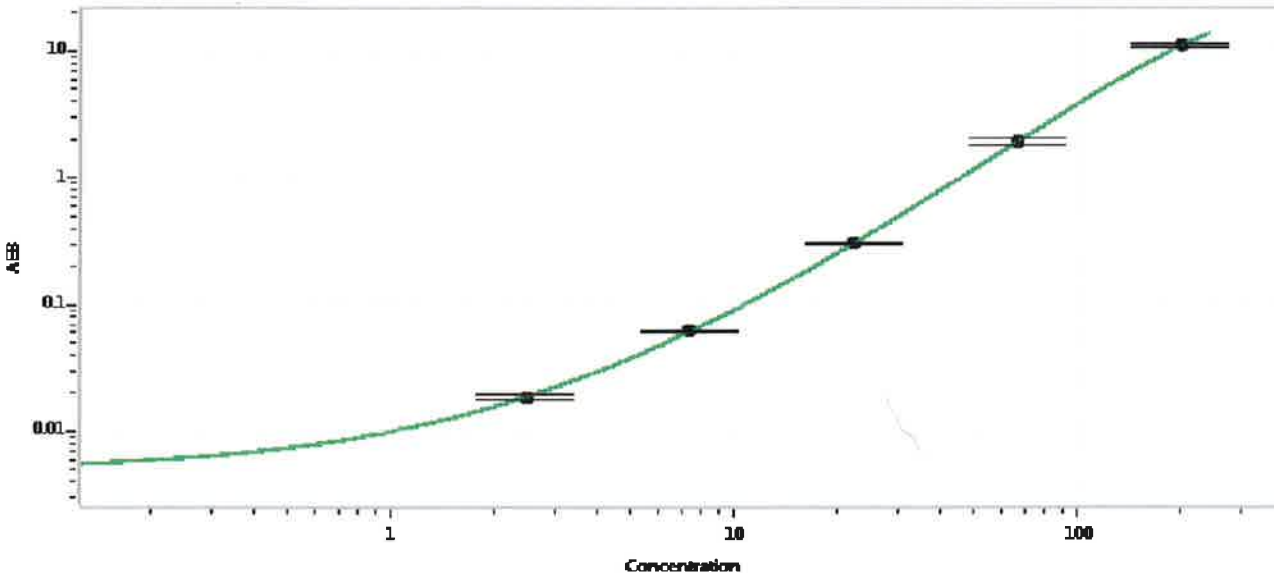
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-021816-02	21.3	17.1-28.3
Control 2	MF-021816-02	321	263-434
Panel 1	MF-012716-05	145	134-226
Panel 2	MF-081415-01	27.8	24.0-39.5

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

Calibrator Concentrate Value Assignment (pg/mL)	2711
---	------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	2.47
C	7.41
D	22.2
E	66.7
F	200

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]*  
Signature/Date  
10 APR 2018