

Product Number:	101599
Lot Number:	501121
Expiration:	9-Aug-2018
Platform(s):	HD-1 & SR-X



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

Component	Item Number	Lot Number
Bead Reagent	101601	73451
Detector Reagent	102012	73452
SBG Reagent	101602	73453
Sample Diluent	101603	73450
Calibrator Diluent	102014	73332
Calibrator Concentrate	100778	73457
RGP Reagent	103159	N/A <sup>1</sup>

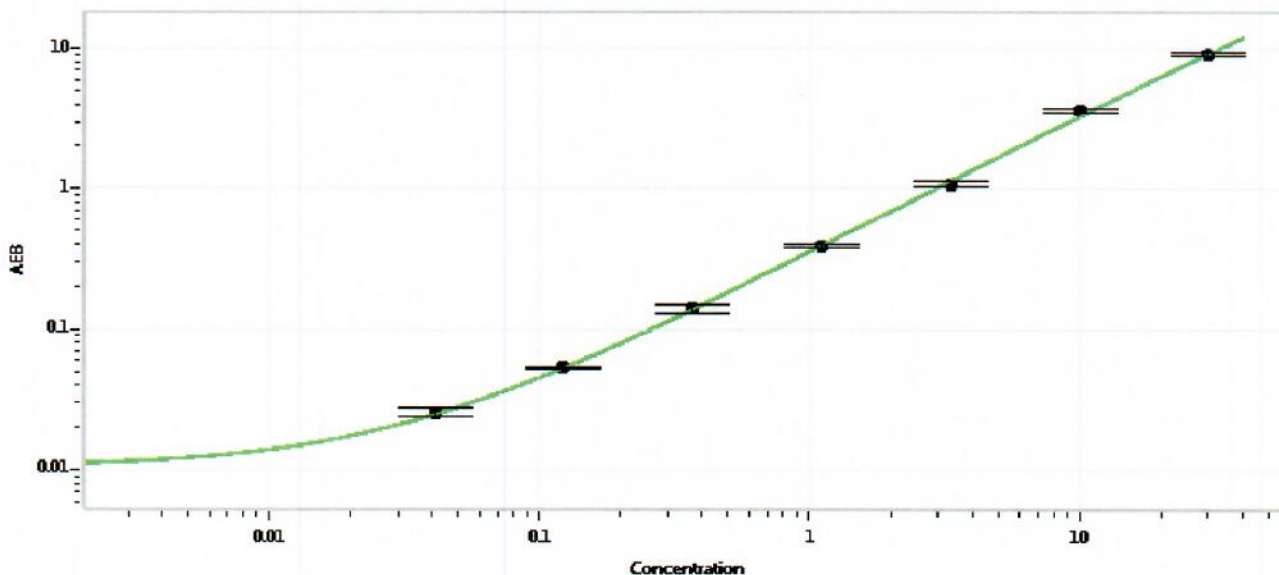
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	52442	0.642	0.485-0.727
Control 2	52442	29.9	19.5-34.7
Panel 1	MF-071615-01	0.350	0.325-0.513
Panel 2	MF-071615-02	7.11	5.58-8.37
Panel 3	MF-071615-03	2.02	1.68-2.52

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

Calibrator Concentrate Value Assignment (pg/mL)	852
---	-----

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

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]*  
Signature/Date  
08 FEB 2018