

Product Number:	101599
Lot Number:	501155
Expiration:	1-Jun-2018
Platform(s):	HD-1 & SR-X



Component	Item Number	Lot Number
Bead Reagent	101601	72492
Detector Reagent	102012	72493
SBG Reagent	101602	72494
Sample Diluent	101603	72441
Calibrator Diluent	102014	72443
Calibrator Concentrate	100778	70764
RGP Reagent	103159	N/A ¹

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

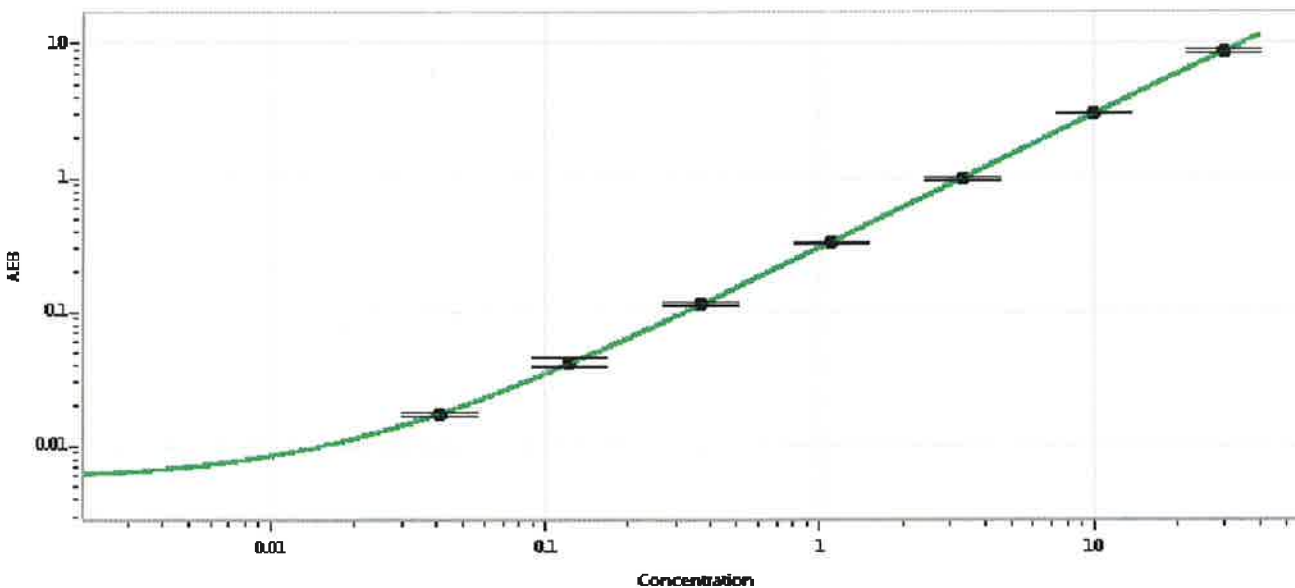
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	52442	0.633	0.485-0.727
Control 2	52442	29.3	19.5-34.7
Panel 1	MF-071615-01	0.440	0.325-0.513
Panel 2	MF-071615-02	8.14	5.58-8.37
Panel 3	MF-071615-03	2.38	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)	636
---	-----

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] 27 FEB 2018
Signature/Date