

Product Number:	101664
Lot Number:	501190
Expiration:	13-Nov-2018
Platform(s):	HD-1



Component	Part Number	Lot Number
Bead Reagent	101666	73382
Detector Reagent	101665	73170
SBG Reagent	101667	73353
Sample Diluent	101668	73171
Calibrator Diluent	101669	73172
Calibrator Concentrate	100717	71660
RGP Reagent	103159	N/A ¹

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

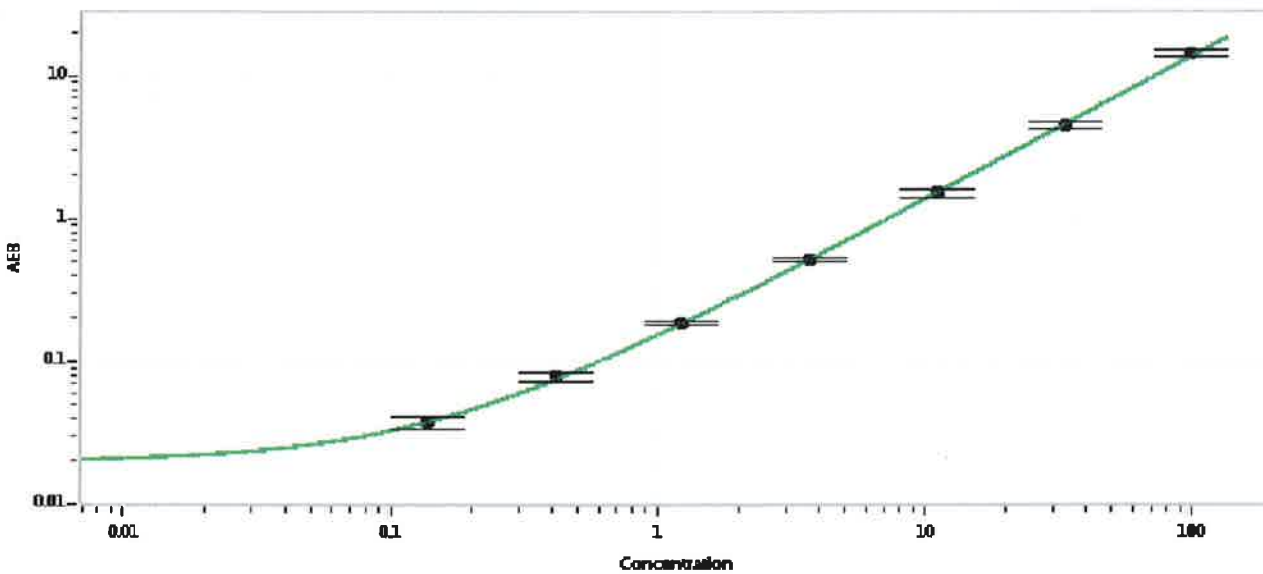
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	52520	1.16	0.740-1.33
Control 2	52520	38.7	27.1-40.6
Panel 1	MF-082015-01	1.06	0.942-1.53
Panel 2	MF-081415-02	9.35	7.46-11.2
Panel 3	MF-081415-03	10.3	9.20-15.0
Panel 4	MF-090315-09	47.2	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)	1991
---	------

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

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 29 MAR 2018
Signature/Date