

Product Number:	101664
Lot Number:	501260
Expiration:	20-Apr-2019
Platform(s):	HD-1 & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101666	81100
Detector Reagent	101665	81136
SBG Reagent	101667	81132
Sample Diluent	101668	81131
Calibrator Diluent	101669	81130
Calibrator Concentrate	100717	71660
RGP Reagent	103159	N/A ¹

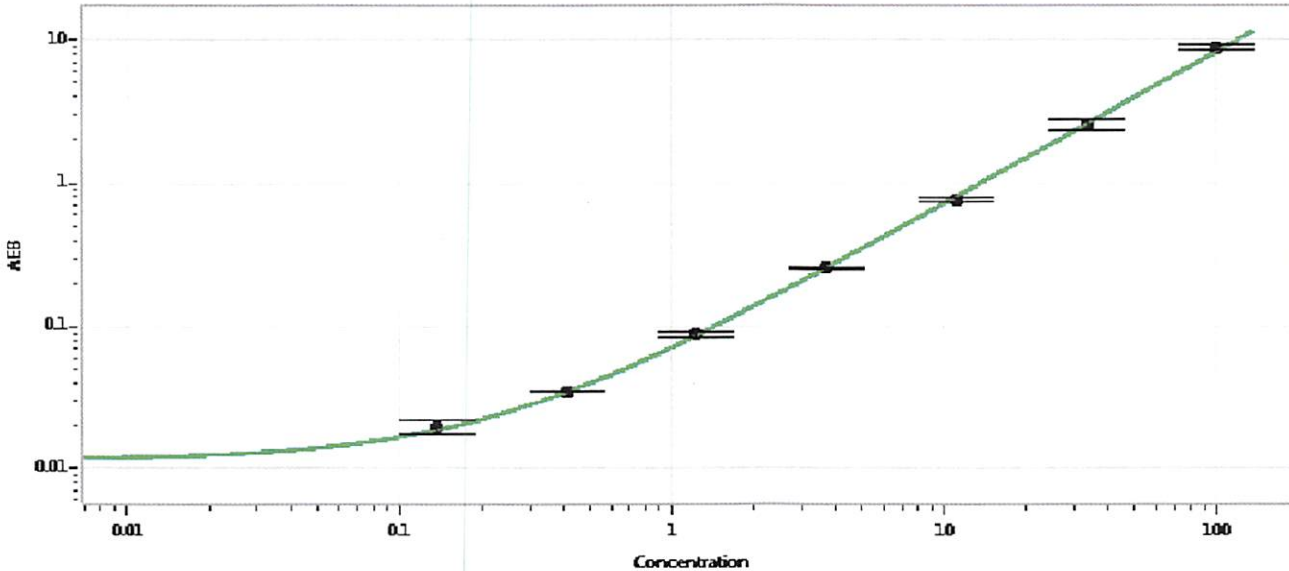
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 2	52520	35.6	27.1-40.6
Panel 1	MF-082015-01	1.30	0.942-1.53
Panel 2	MF-081415-02	8.98	7.46-11.2
Panel 3	MF-081415-03	10.9	9.20-15.0
Panel 4	MF-090315-09	45.4	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] 23 MAY 2018
Signature/Date