

Product Number:	102258
Lot Number:	501102
Expiration:	15-Sep-2018
Platform(s):	HD-1 & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	102246	80151
Detector Reagent	102248	73497
SBG Reagent	102250	73491
Sample Diluent	102252	73495
Calibrator Diluent	102620	73492
Calibrator Concentrate	102619	73520
RGP Reagent	103159	N/A ¹

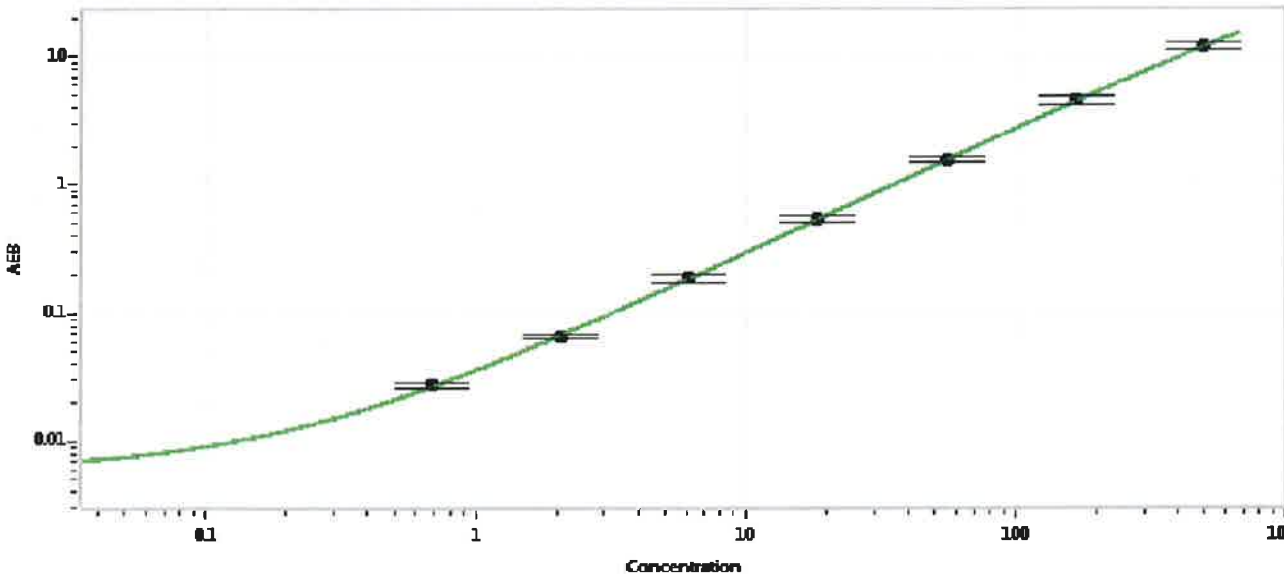
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	72720	5.32	3.87-5.80
Control 2	72720	195	114-225
Panel 1	MF-092717-01	5.25	4.79-7.19
Panel 2	MF-092717-02	59.3	49.7-77.6
Panel 3	MF-092717-03	232	122-297

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

Calibrator Concentrate Value Assignment (pg/mL)	7,585
---	-------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.686
C	2.06
D	6.17
E	18.5
F	55.6
G	167
H	500

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date
06 FEB 2018