

Product Number:	101622
Lot Number:	501376
Expiration:	10-Jan-2019
Platform(s):	HD-1 & SR-X



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

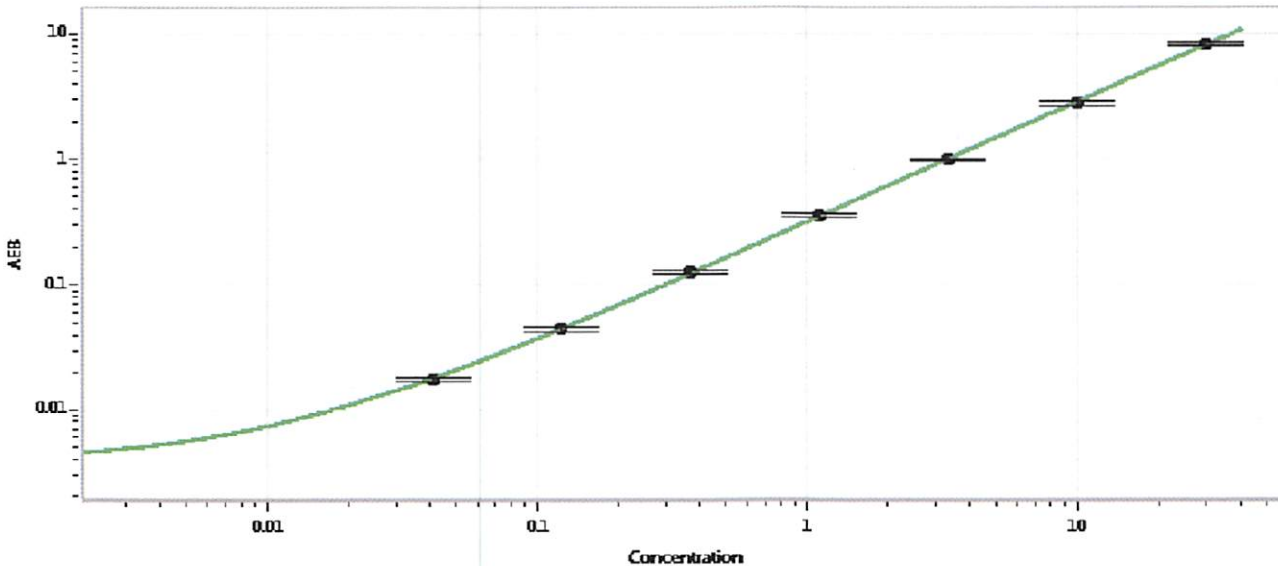
Component	Part Number	Lot Number
Bead Reagent	101624	821504
Detector Reagent	101623	81712
SBG Reagent	101625	822801
Sample Diluent	103203	81004
Calibrator Diluent	101627	81632
Calibrator Concentrate	100652	72350
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	52173	1.74	1.27-1.91	
Control 2	52173	31.3	19.1-31.4	
Panel 1	MF-050918-01	54.2	40.7-65.6	
Panel 2	MF-050918-02	7.28	5.20-8.12	
Panel 3	MF-050918-03	0.971	0.620-1.21	

Calibrator Concentrate Value Assignment (pg/mL)	797
---	-----

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] 12/19/2018
Signature/Date