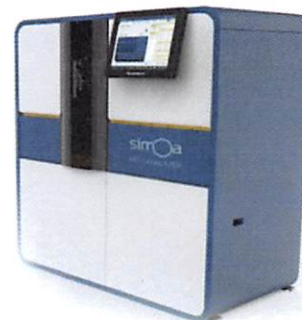


Product Number:	101478
Lot Number:	501433
Expiration:	16-Feb-2019
Platform(s):	HD-1 & SR-X



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

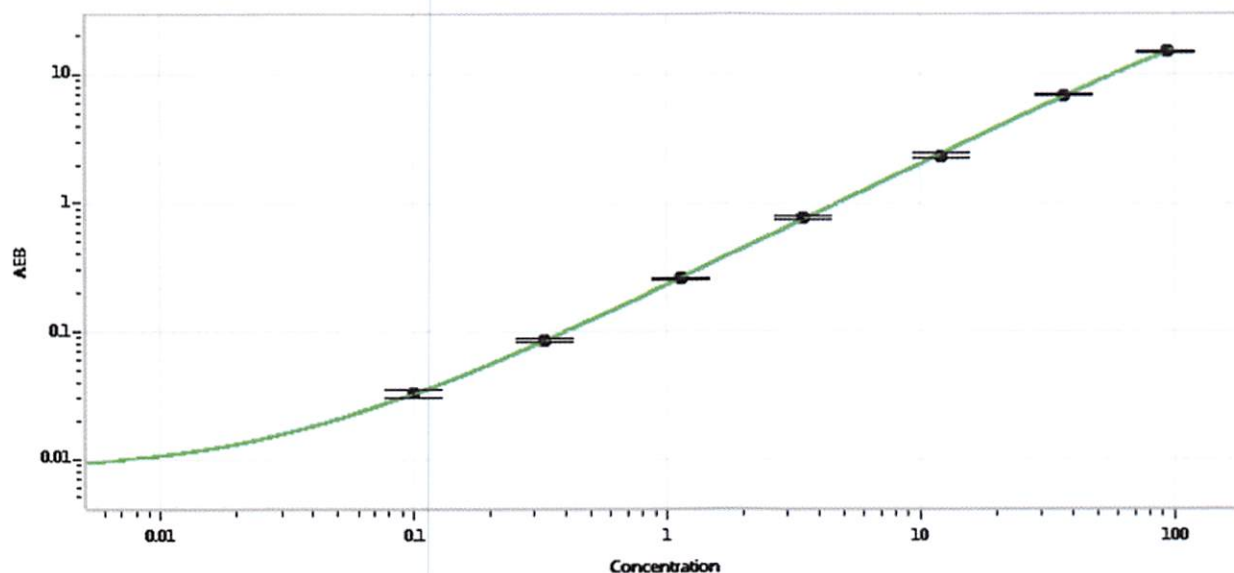
Component	Part Number	Lot Number
Bead Reagent	101480	826705
Detector Reagent	101481	826212
SBG Reagent	101482	822810
Sample Diluent	101483	827001
Calibrators	100758-100765	80230
Controls	100768 & 100769	80231
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80231	3.32	2.56-3.84
Control 2	80231	75.6	58.1-87.2
Panel 1	MF-051017-04	0.732	0.508-0.855
Panel 2	MF-042117-06	5.07	3.89-6.07
Panel 3	MF-042117-07	63.0	45.5-73.7

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.100
C	0.327
D	1.14
E	3.44
F	12.1
G	37.1
H	93.4

Review/Approval

Sivasankari Murugan

QC Supervisor

Name

Title

Signature/Date

*[Handwritten Signature]* 05 OCT 2018