

Product Number:	101478
Lot Number:	501365
Expiration:	3-Feb-2019
Platform(s):	HD-1 & SR-X



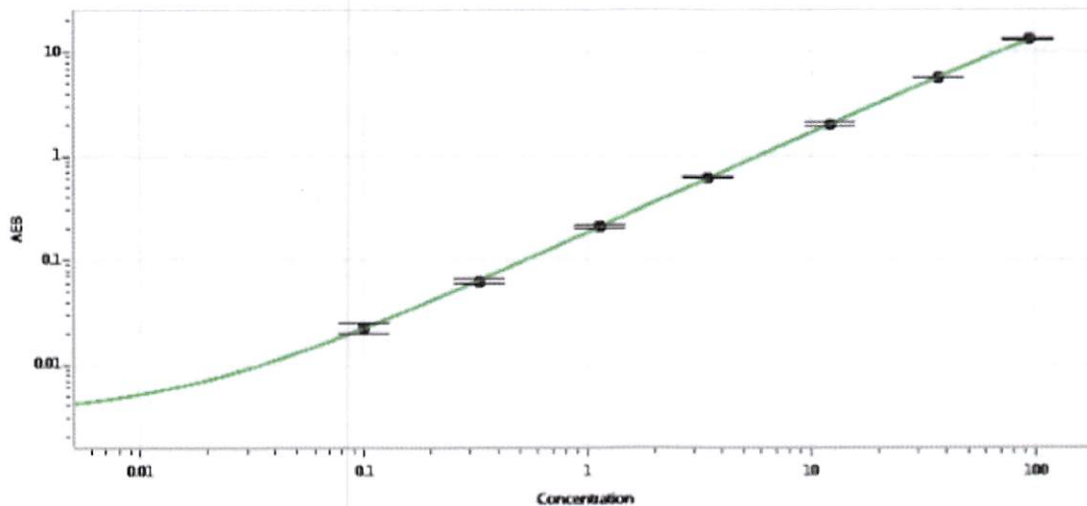
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101480	822001
Detector Reagent	101481	821802
SBG Reagent	101482	821502
Sample Diluent	101483	821907
Calibrators	100758-100765	80230
Controls	100768 & 100769	80231
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	80231	3.27	2.56-3.84	
Control 2	80231	70.0	58.1-87.2	
Panel 1	MF-051017-04	0.660	0.508-0.855	
Panel 2	MF-042117-06	4.40	3.89-6.07	
Panel 3	MF-042117-07	49.3	45.5-73.7	

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] 14 Aug 2018