

Product Number:	103277
Lot Number:	501523
Expiration:	20-May-2019
Platform(s):	HD-1



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

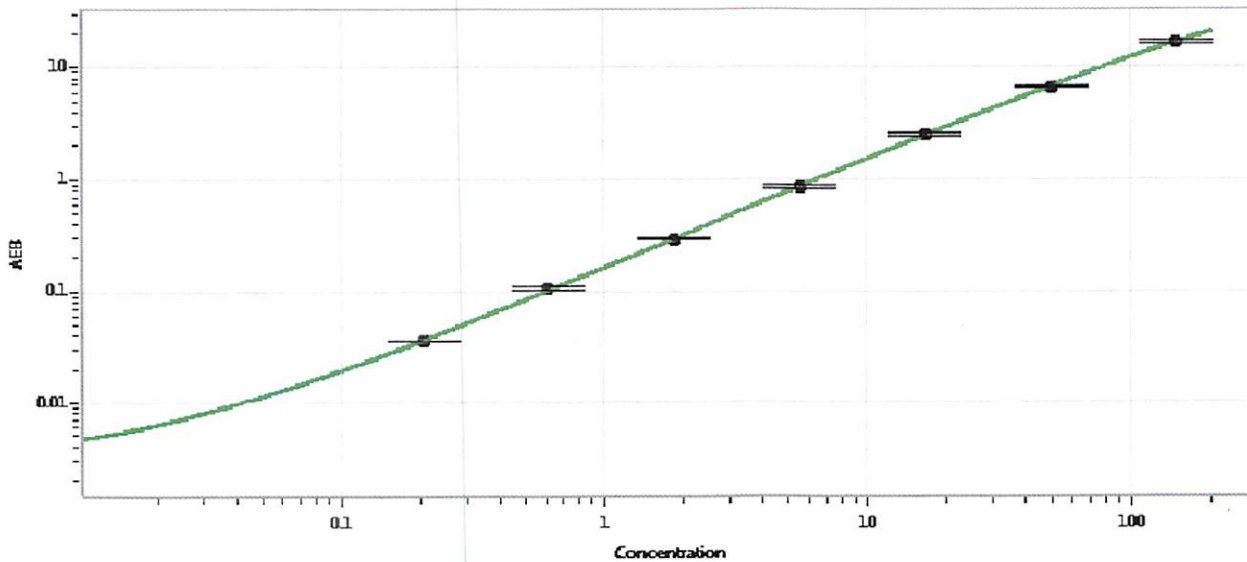
Component	Part Number	Lot Number
Bead Reagent	103265	832404
Detector Reagent	103268	832505
SBG Reagent	103269	830504
Sample Diluent	103270	832402
Calibrator Diluent	103266	830403
Calibrator Concentrate	103272	832501
RGP Reagent	103159	N/A ¹

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	824302	5.01	4.22-6.33	
Control 2	824302	131	111-167	
Panel 1	823501	2.44	1.85-3.04	
Panel 2	823502	30.4	21.3-37.8	
Panel 3	823503	92.9	71.6-107	

Calibrator Concentrate Value Assignment (pg/mL)	5550
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.206
C	0.617
D	1.85
E	5.56
F	16.7
G	50.0
H	150

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date 10 DEC 2018