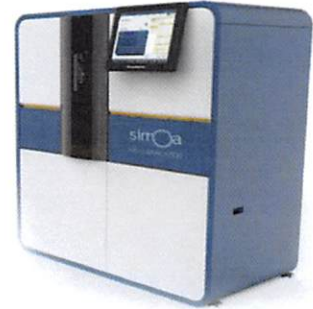


Product Number:	102732
Lot Number:	501462
Expiration:	9-Apr-2019
Platform(s):	HD-1 & SR-X



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

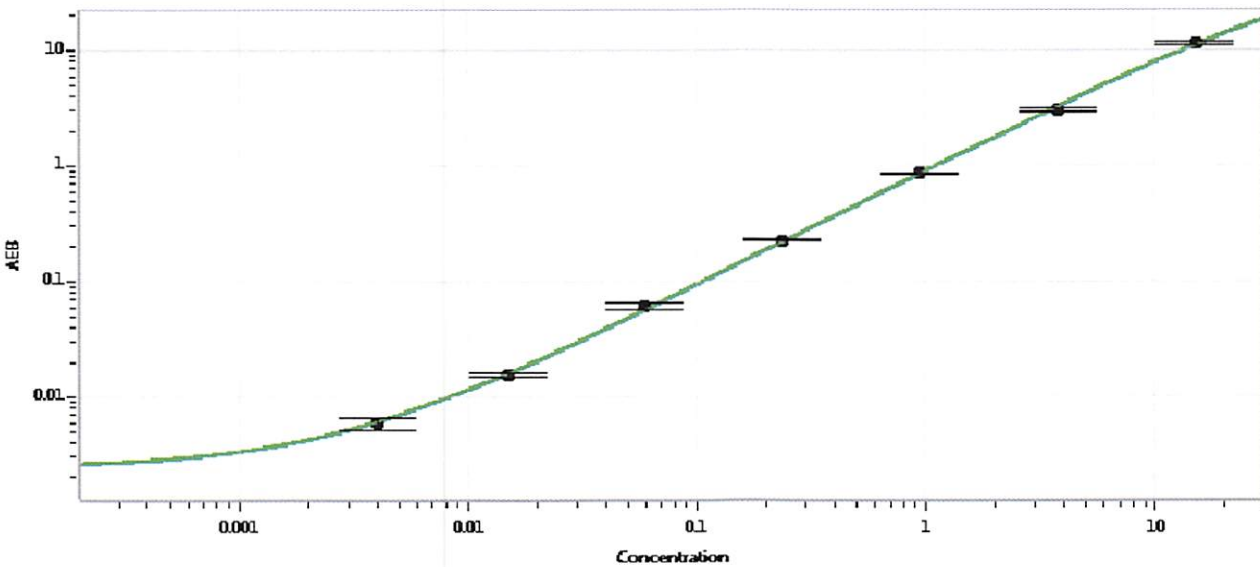
Component	Part Number	Lot Number
Bead Reagent	102722	829007
Detector Reagent	102724	828205
SBG Reagent	102726	828501
Sample Diluent	102728	828503
Calibrator Diluent	102731	826810
Calibrator Concentrate	100658	73217
RGP Reagent	103159	N/A ¹

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	MF-061217-01	0.180	0.134-0.289	
Control 2	MF-061217-01	3.07	2.44-6.35	
Panel 1	MF-061317-03	0.040	0.037-0.088	
Panel 2	MF-061317-04	0.258	0.191-0.490	
Panel 3	MF-061317-05	4.06	3.18-7.89	

Calibrator Concentrate Value Assignment (pg/mL)	194
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.004
C	0.015
D	0.059
E	0.234
F	0.938
G	3.75
H	15.0

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date