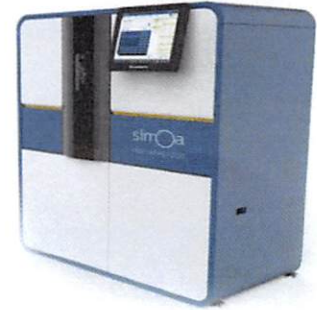


Product Number:	101154
Lot Number:	501435
Expiration:	24-Mar-2019
Platform(s):	HD-1

Component	Part Number	Lot Number
Bead Reagent	101141	827110
Detector Reagent	101143	826703
SBG Reagent	101153	826702
Sample Diluent	103320	826211
Calibrator Diluent	101147	823613
Calibrator Concentrate	101145	73524
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific



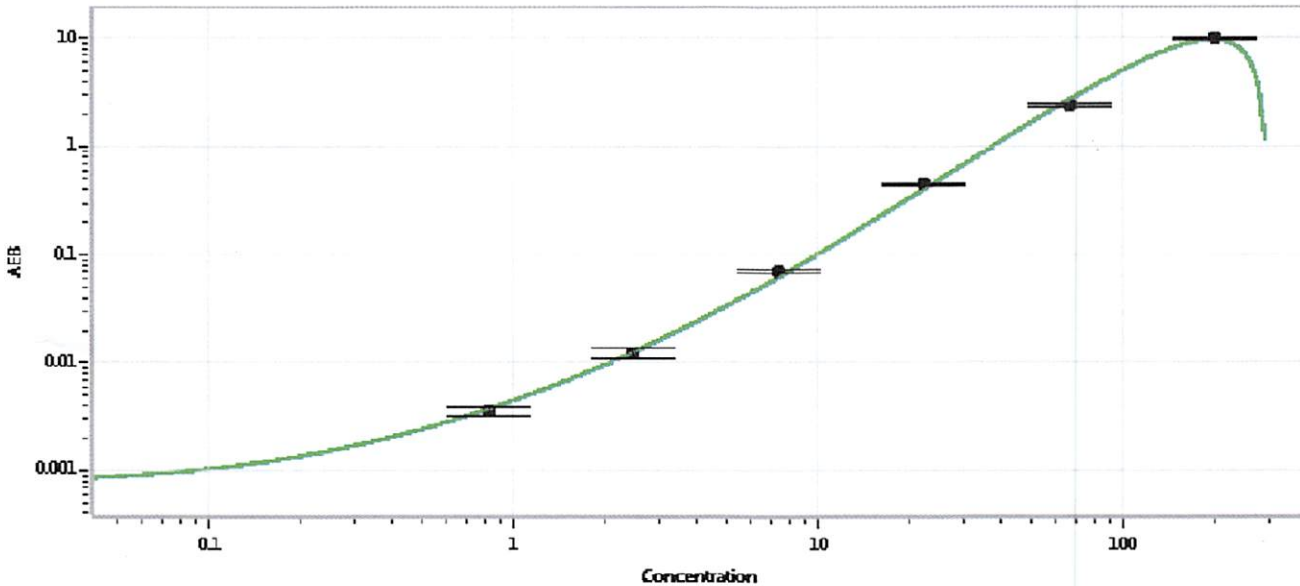
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	41952	68.8	51.8-77.6
Control 2	41952	397	269-404
Panel 1	1816	105	75.5-113
Panel 2	1712	99.9	72.8-109
Panel 3	2004	159	114-170

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

Calibrator Concentrate Value Assignment (pg/mL)	14,199
-------------------------------------------------	--------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.823
C	2.47
D	7.41
E	22.2
F	66.7
G	200

Review/Approval: Sivasankari Murugan QC Supervisor [Signature] 09 OCT 2018
Name Title Signature/Date