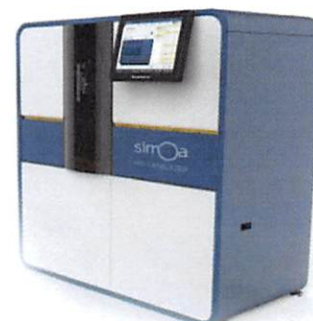


Product Number:	102669
Lot Number:	501423
Expiration:	11-Mar-2019
Platform(s):	HD-1



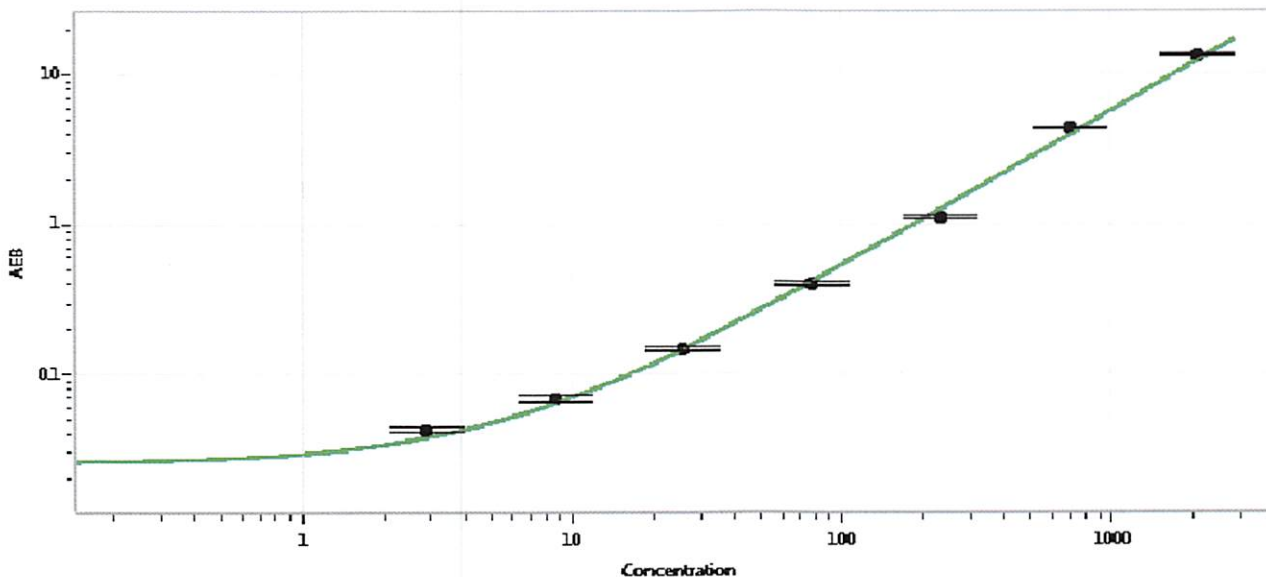
Data below represents results generated on the Simoa™ HD-1

Component	Part Number	Lot Number
Bead Stock	102662	825508
Detector Stock	102663	825509
Calibrator Concentrate	102666	MF-020718-03
Calibrator Diluent	102664	825406
Sample Diluent	102665	825407

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159

Calibrator Concentrate Target Value	60 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.88
C	8.64
D	25.9
E	77.8
F	233
G	700
H	2100

Review/Approval	Sivasankari Murugan	QC Supervisor	<i>B. Murugan</i> 24 Oct 2018
	Name	Title	Signature/Date