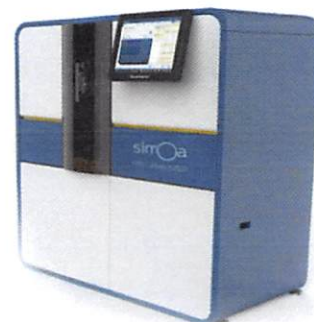


Product Number:	102669
Lot Number:	501790
Expiration:	19-Jan-2020
Platform(s):	HD-1



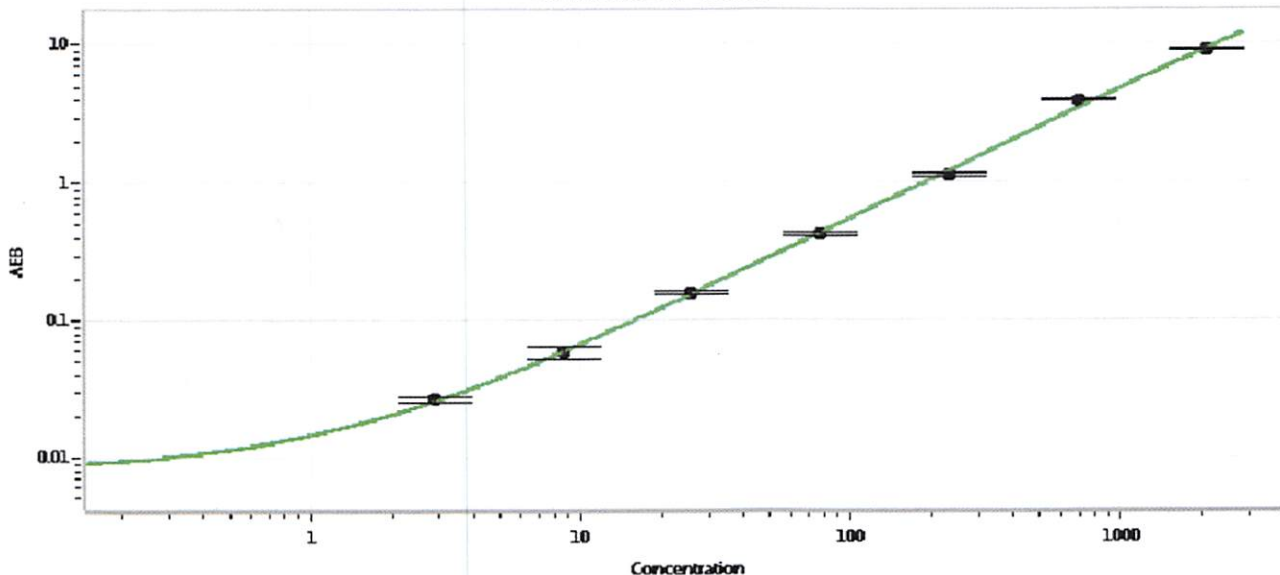
Data below represents results generated on the Simoa™ HD-1

Component	Part Number	Lot Number
Bead Stock	102662	902405
Detector Stock	102663	902505
Calibrator Concentrate	102666	MF-020718-03
Calibrator Diluent	102664	901911
Sample Diluent	102665	902506

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		101957
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

Patrizia Stadler
Name

QC Assoc. Director
Title

p. stadler 26 Apr 2019
Signature/Date