

Product Number:	102491
Lot Number:	501710
Expiration:	7-Sep-2019
Platform(s):	HD-1

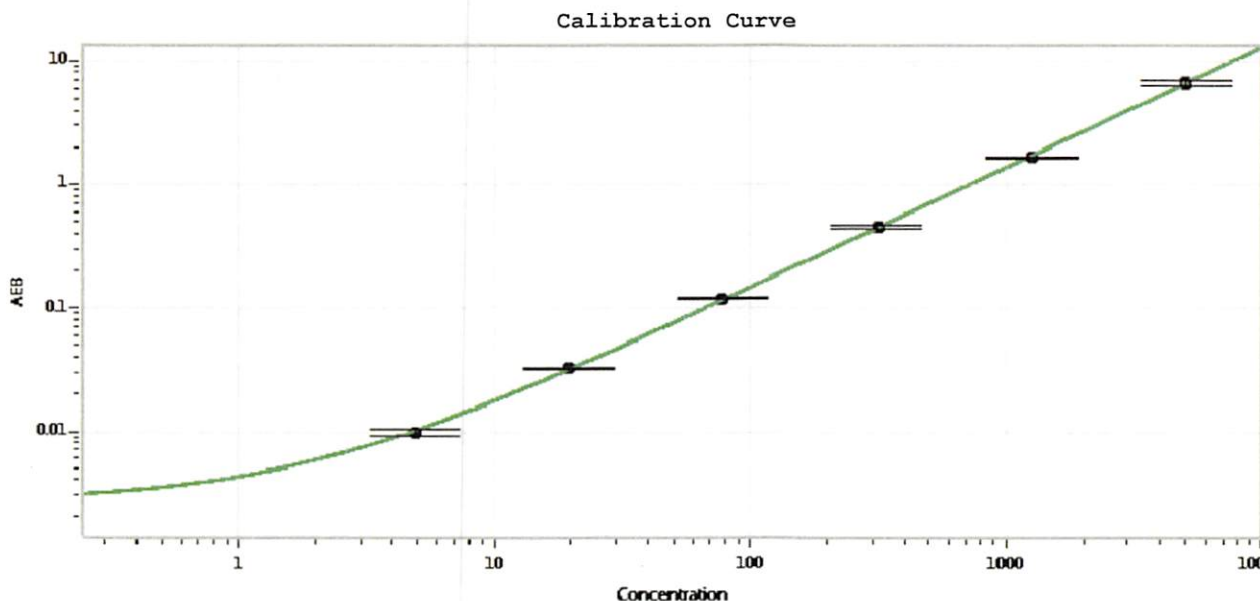


Data below represents results generated on the Simoa™ HD-1

Component	Part Number	Lot Number
Bead Stock	102484	906705
Detector Stock	102485	906706
Detector Diluent	103084	906613
Calibrator Concentrate	102488	907004
Calibrator Diluent	102486	906612
Sample Diluent	102487	906611

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

Calibrator Concentrate Target Value	100 ng/mL
-------------------------------------	-----------



Calibrator Levels (pg/mL)	
A	0.000
B	4.88
C	19.5
D	78.1
E	313
F	1250
G	5000

Review/Approval

Patrizia Stadler  
Name

Assoc. Director, QC  
Title

*P. Stadler* 14 Mar 2019  
Signature/Date