

Product Number:	101419
Lot Number:	501775
Expiration:	21-Dec-19
Platform(s):	HD-1



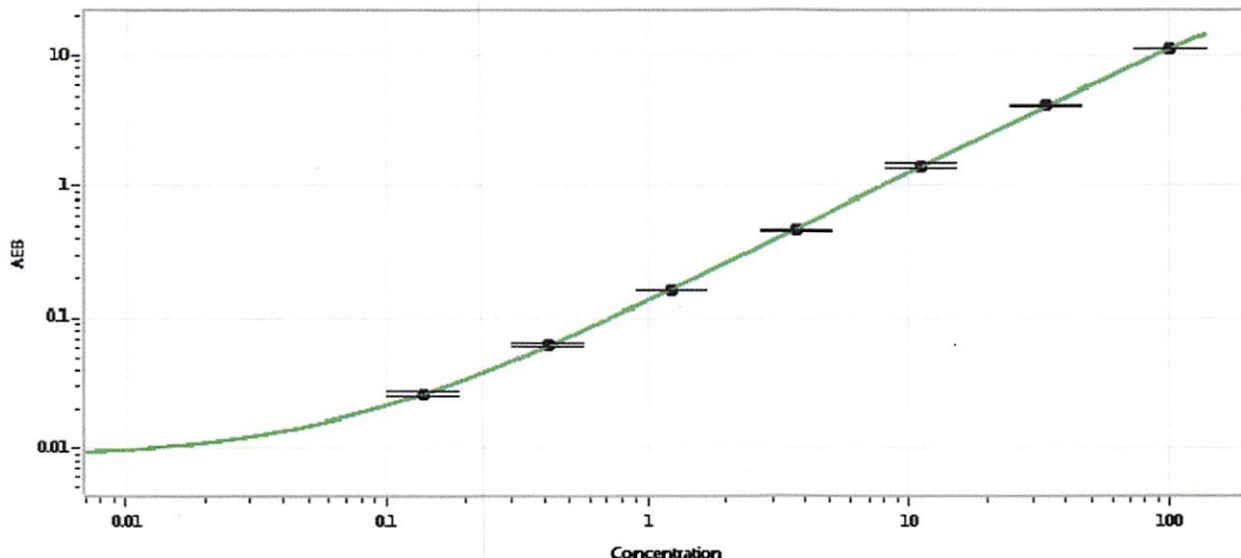
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	101899	910808
Detector Stock	101900	910502
Calibrator Concentrate	101420	828601
Calibrator Diluent	101418	908820
Sample Diluent	101416	835509

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

Calibrator Concentrate Target Value	2 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.137
C	0.412
D	1.23
E	3.70
F	11.1
G	33.3
H	100

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	24 Apr 2019
	Name	Title	Signature/Date	