

Product Number:	102343
Lot Number:	501725
Expiration:	21-Dec-2019
Platform(s):	HD-1



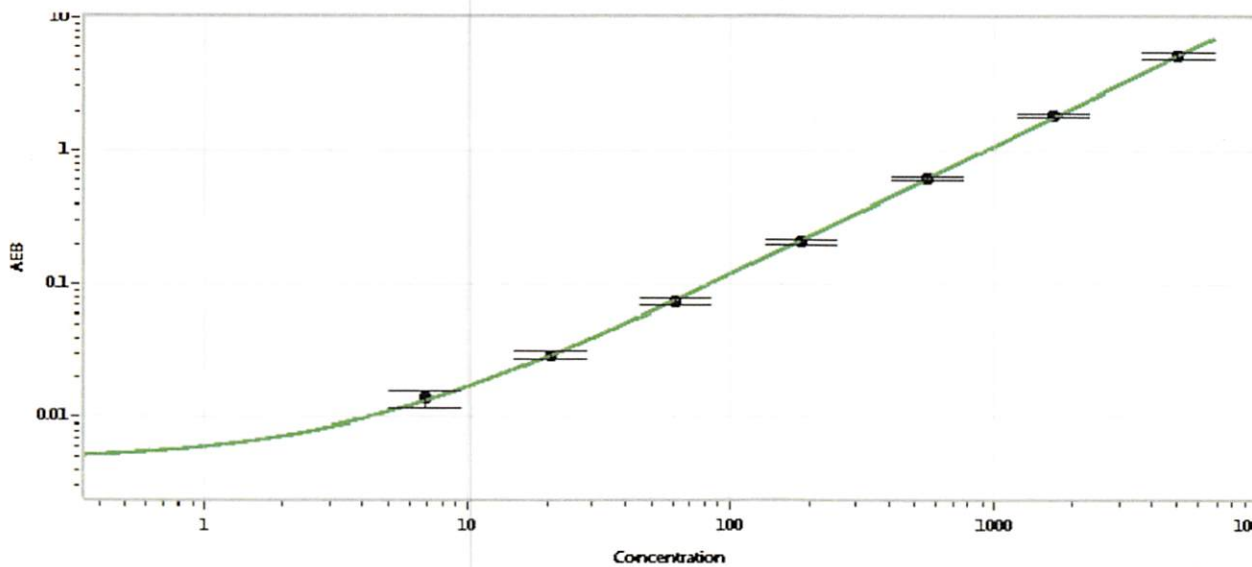
Data below represents results generated on the Simoa™ HD-1

Component	Part Number	Lot Number
Bead Stock	102337	907707
Detector Stock	102338	907301
Calibrator Concentrate	102342	831815
Calibrator Diluent	102339	835506
Sample Diluent	102340	907008

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	100 ng/mL
-------------------------------------	-----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	6.86
C	20.6
D	61.7
E	185
F	556
G	1667
H	5000

Review/Approval: Patrizia Stadler Assoc. Director QC P. Stadler 28 Mar 2019
 Name Title Signature/Date