

Product Number:	101871
Lot Number:	502033
Expiration:	30-Mar-2020
Platform(s):	HD-1/HD-X

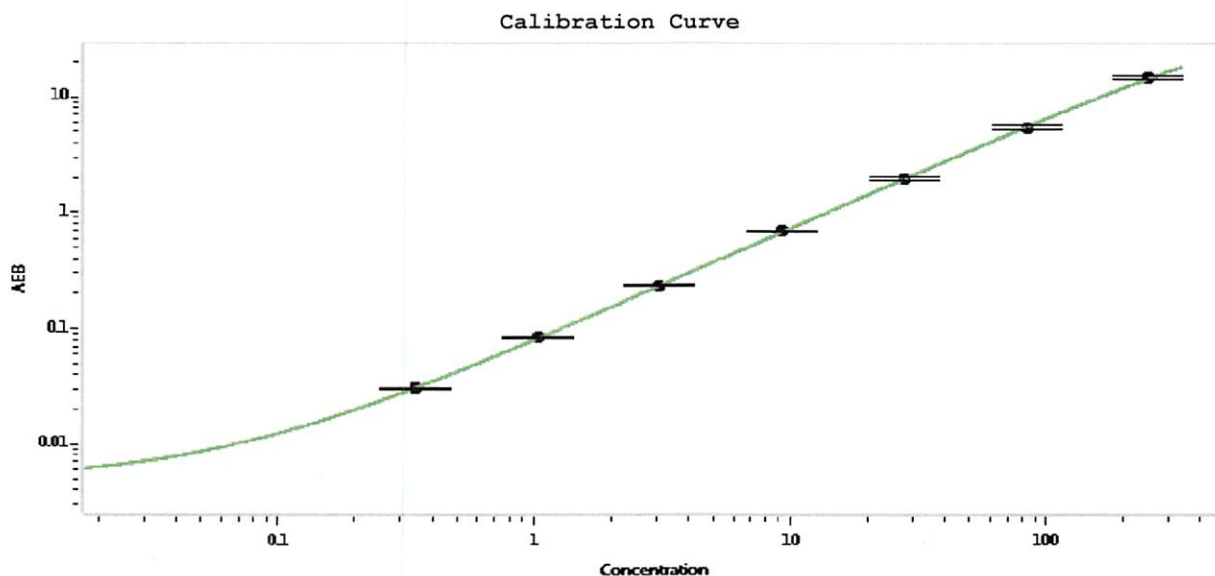


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	101893	926012
Detector Stock	101894	926008
Calibrator Concentrate	101866	926009
Calibrator Diluent	101959	924904
Sample Diluent	101958	908920

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	5 ng/mL
-------------------------------------	---------



Calibrator Levels (pg/mL)	
A	0.000
B	0.343
C	1.03
D	3.09
E	9.26
F	27.8
G	83.3
H	250

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	25 Sep 2019
	Name	Title	Signature/Date	