

Product Number:	102336
Lot Number:	502028
Expiration:	30-Mar-2020
Platform(s):	HD-1/HD-X



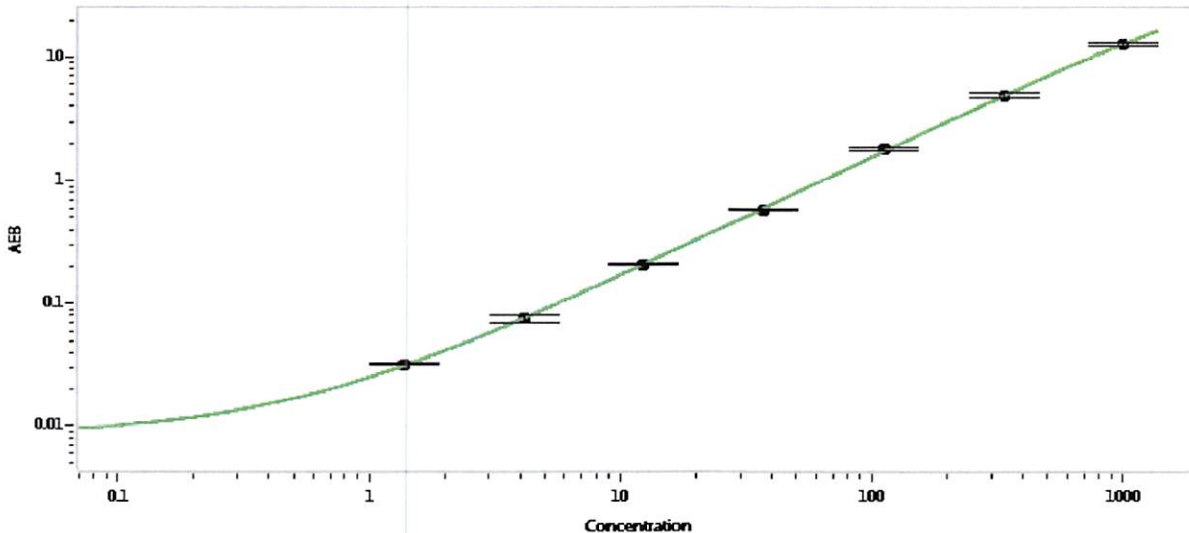
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102330	925208
Detector Stock	102331	925507
Calibrator Concentrate	102335	920503
Calibrator Diluent	102332	908918
Sample Diluent	102333	925407

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	20 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.37
C	4.11
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	18 Sep 2019
	Name	Title	Signature/Date	