

Product Number:	102583
Lot Number:	501783
Expiration:	30-Mar-2020
Platform(s):	HD-1

Component	Part Number	Lot Number
Bead Stock	102576	911206
Detector Stock	102577	911409
Calibrator Concentrate	102580	912804
Calibrator Diluent	102578	910307
Sample Diluent	102579	908904

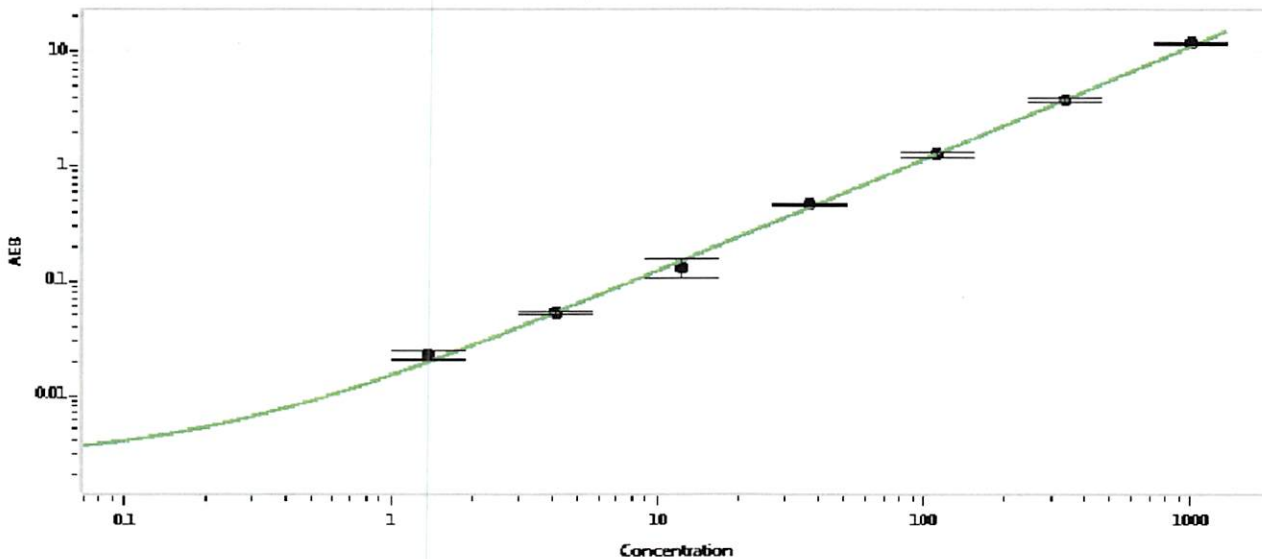


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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	10 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.37
C	4.12
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval	Patrizia Stadler	QC Asso. Director	<i>P. Stadler</i>	16 July 2019
	Name	Title	Signature/Date	