

Product Number:	102583
Lot Number:	501666
Expiration:	16-May-2019
Platform(s):	HD-1

Component	Part Number	Lot Number
Bead Stock	102576	820810
Detector Stock	102577	819710
Calibrator Concentrate	102580	MF-051617-03
Calibrator Diluent	102578	816309
Sample Diluent	102579	819705

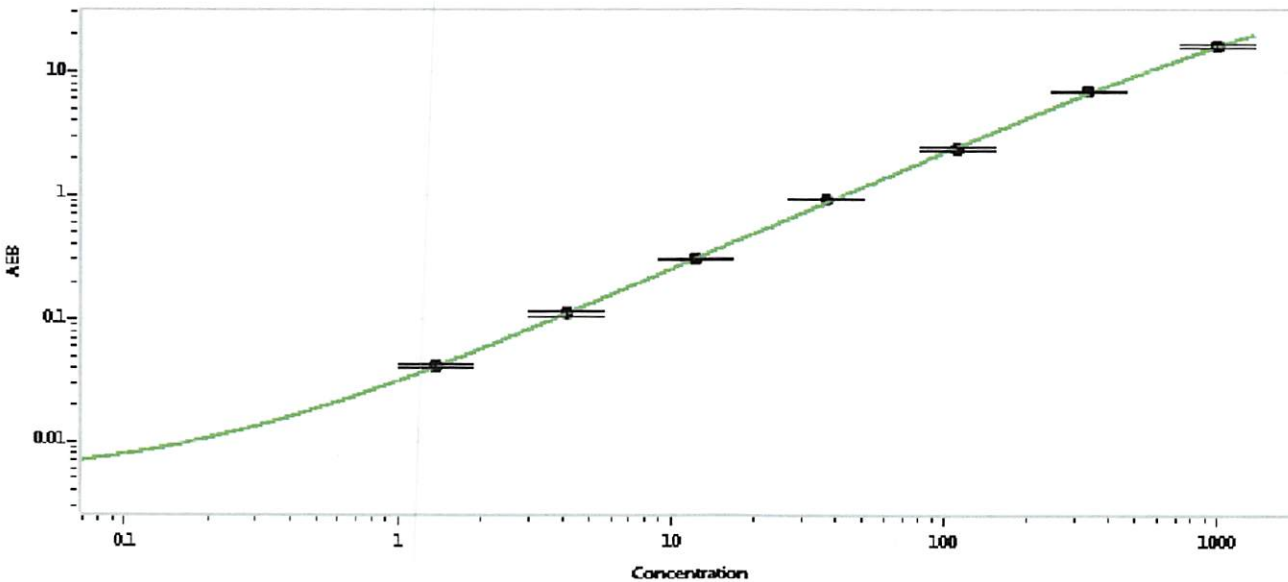


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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>	15 Feb 2019
	Name	Title	Signature/Date	