

Product Number:	103093
Lot Number:	501916
Expiration:	8-Jan-2020
Platform(s):	HD-1/ HD-X & SR-X



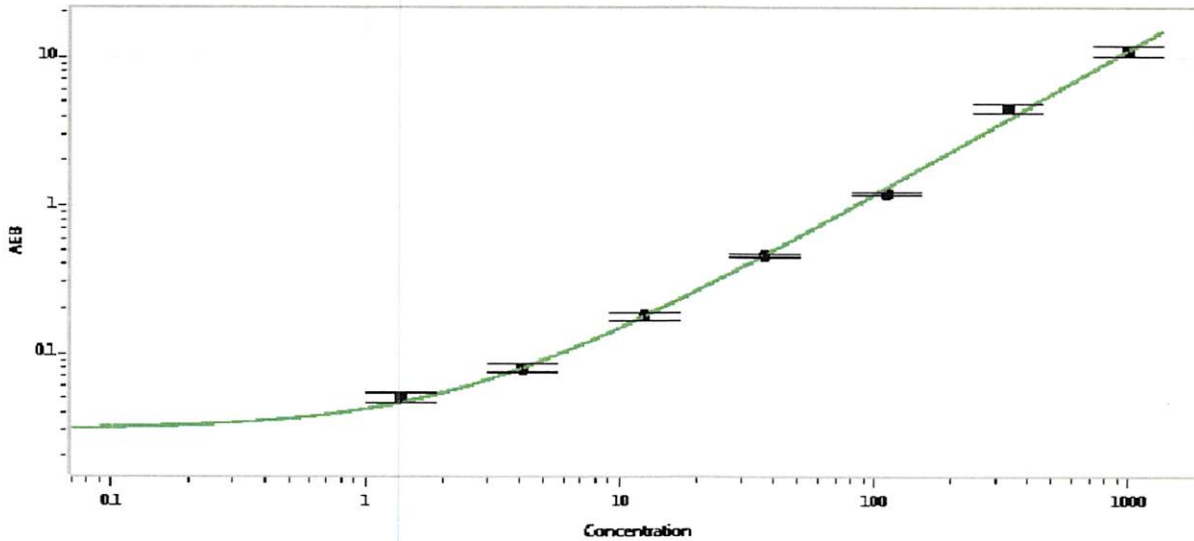
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102021	912206
Detector Stock	103043	919301
Calibrator Concentrate	103047	MF-010818-10
Calibrator Diluent	103044	912203
Sample Diluent	103046	912202

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	12 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	Assoc. Director, QC	<i>P. Stadler</i> 17 Jul 2019
	Name	Title	Signature/Date