

Product Number:	103093
Lot Number:	502196
Expiration:	2-May-2020
Platform(s):	HD-1/ HD-X & SR-X

Component	Part Number	Lot Number
Bead Stock	102021	912206
Detector Stock	103043	919301
Calibrator Concentrate	103047	935308
Calibrator Diluent	103044	912203
Sample Diluent	103046	912202

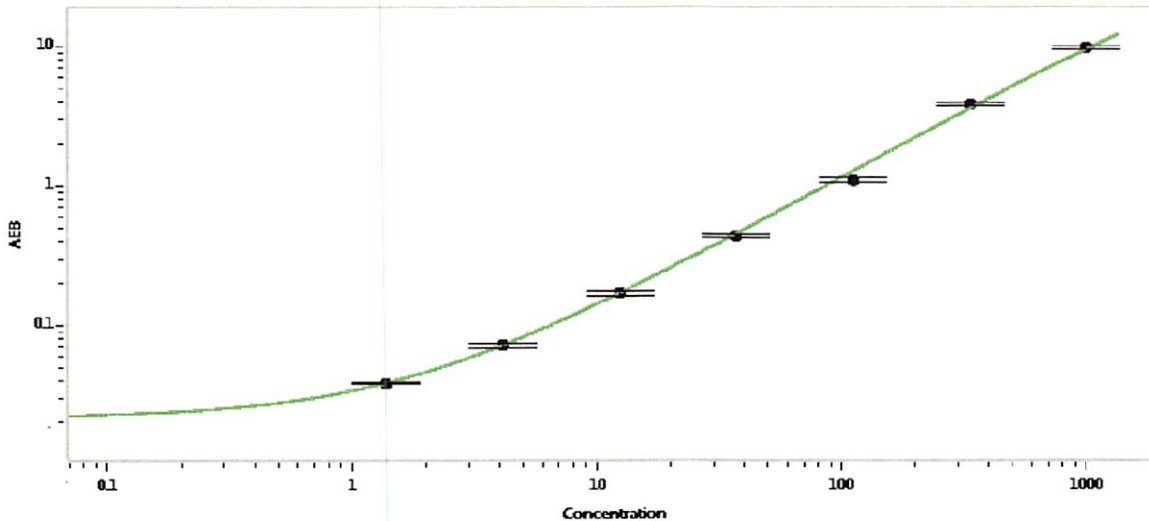


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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	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  
Name

Assoc. Director, QC  
Title

*p.stadler 2/10/20*  
Signature/Date