

Product Number:	103071
Lot Number:	501740
Expiration:	29-Jan-2020
Platform(s):	HD-1 & SR-X

Component	Part Number	Lot Number
Bead Stock	103064	908106
Detector Stock	103065	908102
Calibrator Concentrate	103068	MF-020418-01
Calibrator Diluent	103066	905823
Sample Diluent	103067	902935



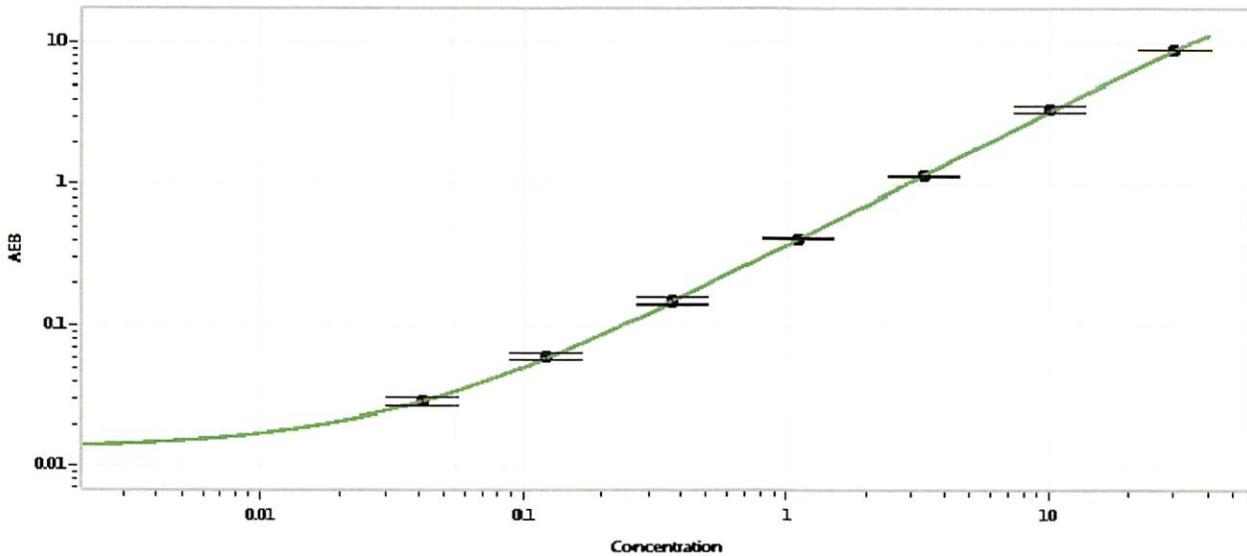
Data below represents results generated on the Simoa™ HD-1 Analyzer.

**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	600 pg/mL
-------------------------------------	-----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval

Patrizia Stadler Assoc. Director, QC  
Name Title

*P Stadler* 29 Mar 2019  
Signature/Date