

Product Number:	103071
Lot Number:	501685
Expiration:	12-Jun-2019
Platform(s):	HD-1 & SR-X

Component	Part Number	Lot Number
Bead Stock	103064	818421
Detector Stock	103065	818422
Calibrator Concentrate	103068	MF-020418-01
Calibrator Diluent	103066	816311
Sample Diluent	103067	818003

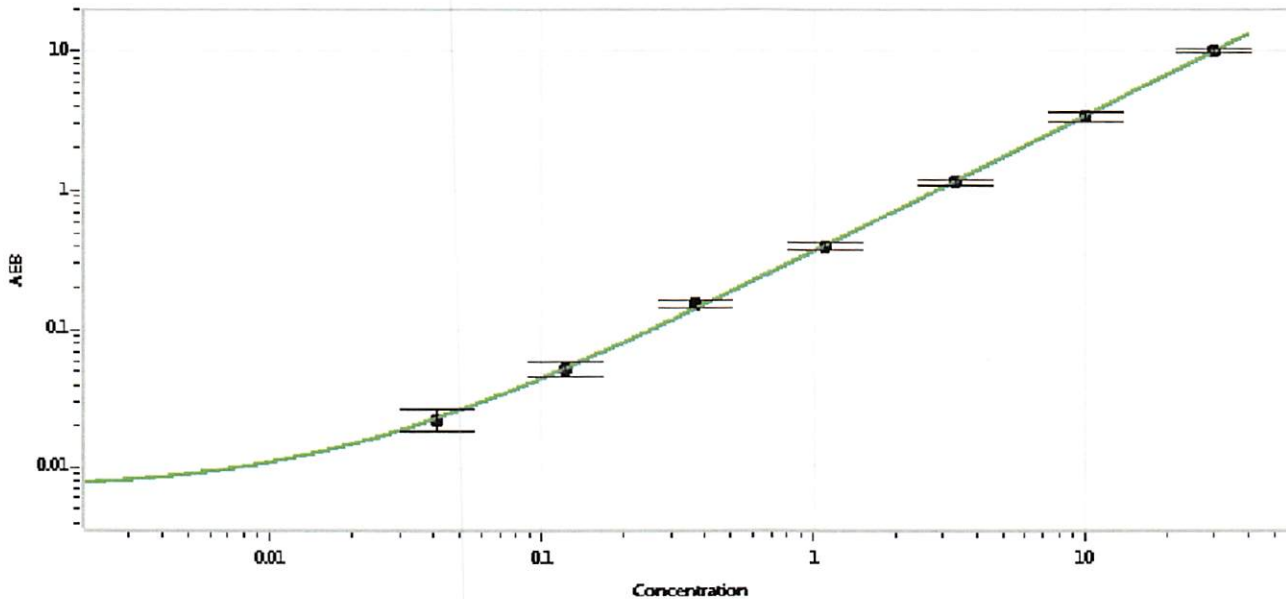


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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 04 Nov 2019
Signature/Date