

Product Number:	103071
Lot Number:	502133
Expiration:	13-May-2020
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Stock	103064	913502
Detector Stock	103065	913309
Calibrator Concentrate	103068	934309
Calibrator Diluent	103066	930820
Sample Diluent	103067	927722

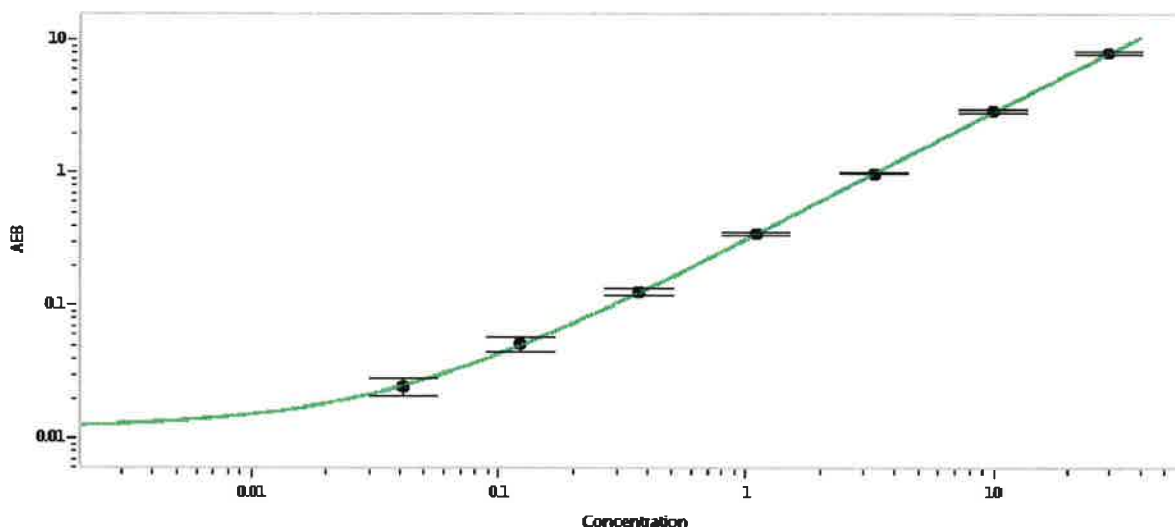


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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
Name

Assoc. Director, QC
Title

P. Stadler 16 Dec 2019
Signature/Date