

Product Number:	102039
Lot Number:	502129
Expiration:	17-Sep-2020
Platform(s):	HD-1/HD-X & SR-X



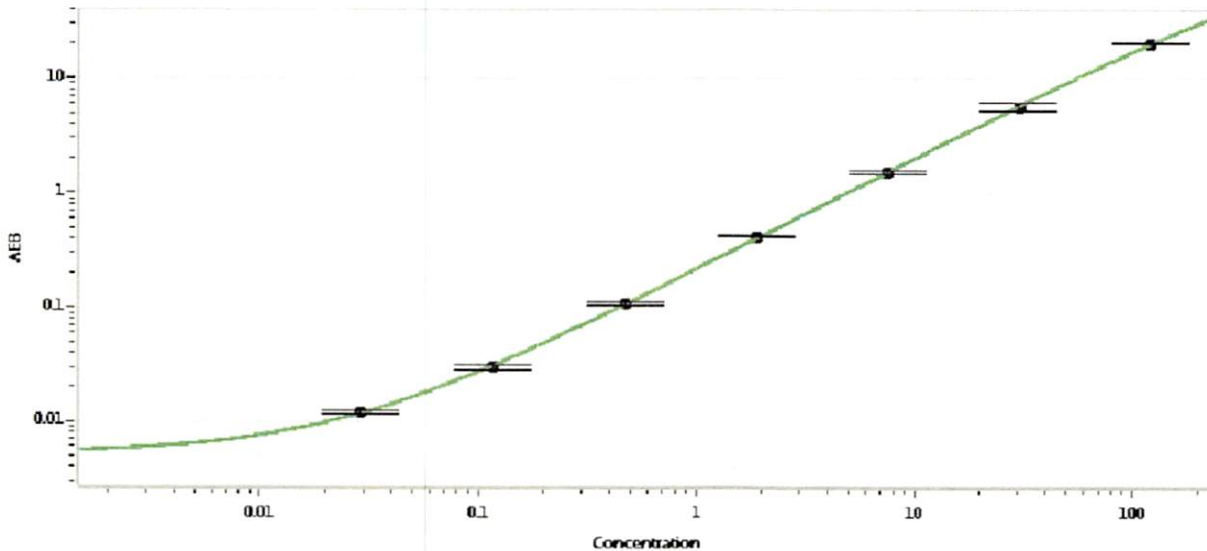
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102028	926205
Detector Stock	102029	926201
Calibrator Concentrate	102032	910804
Calibrator Diluent	102030	926013
Sample Diluent	103402	927719

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	2 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.029
C	0.117
D	0.469
E	1.88
F	7.50
G	30.0
H	120

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

p. Stadler 09 Dec 2019
Signature/Date