

Product Number:	102318
Lot Number:	502079
Expiration:	26-Mar-2020
Platform(s):	HD-1/HD-X

Component	Part Number	Lot Number
Bead Stock	102309	929002
Detector Stock	102310	908520
Calibrator Concentrate	102314	929702
Calibrator Diluent	102311	929102
Sample Diluent	102312	929101

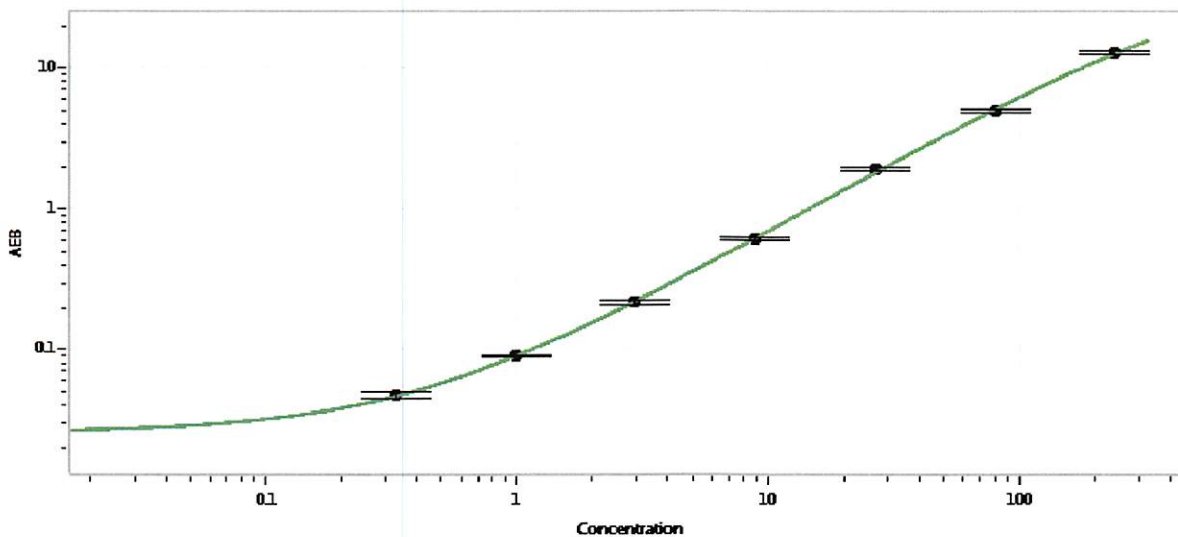


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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	4,800 pg/mL
-------------------------------------	-------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.329
C	0.988
D	2.96
E	8.89
F	26.7
G	80
H	240

Review/Approval

Patrizia Stadler  
Name

Assoc. Director, QC  
Title

*P. Stadler* 5 Nov 2019  
Signature/Date