

Product Number:	102669
Lot Number:	501978
Expiration:	13-May-2020
Platform(s):	HD-1/HD-X



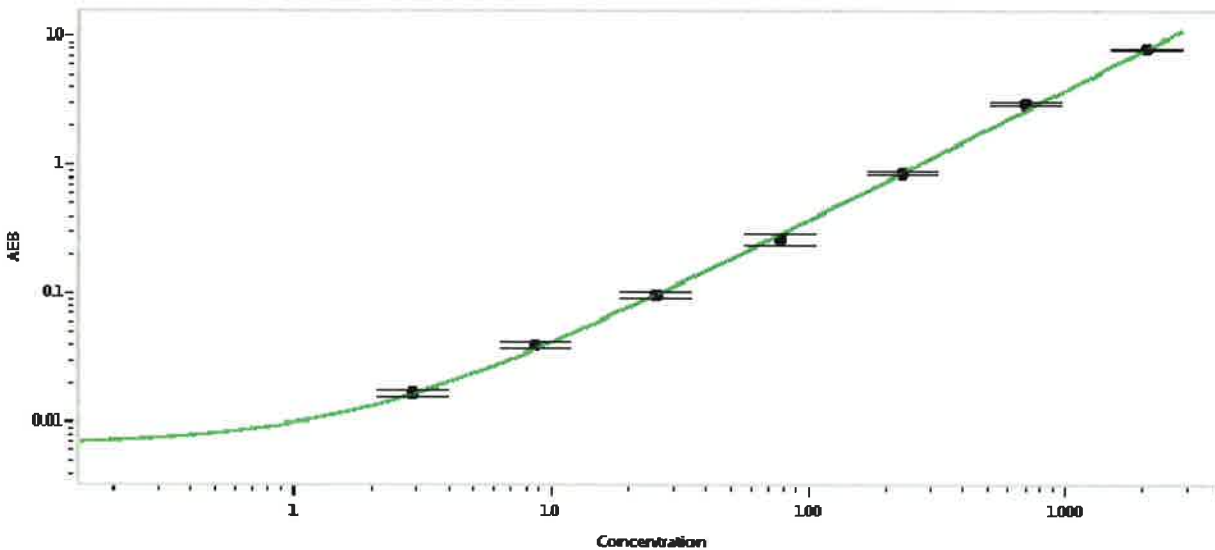
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102662	921023
Detector Stock	102663	913310
Calibrator Concentrate	102666	913401
Calibrator Diluent	102664	922105
Sample Diluent	102665	921312

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	60 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.88
C	8.64
D	25.9
E	77.8
F	233
G	700
H	2100

Review/Approval

Patrizia Stadler  
Name

Assoc. Director, QC  
Title

*P. Stadler* 12 Aug 2019  
Signature/Date