

Product Number:	102209
Lot Number:	501969
Expiration:	19-Jun-2020
Platform(s):	HD-1/HD-X & SR-X



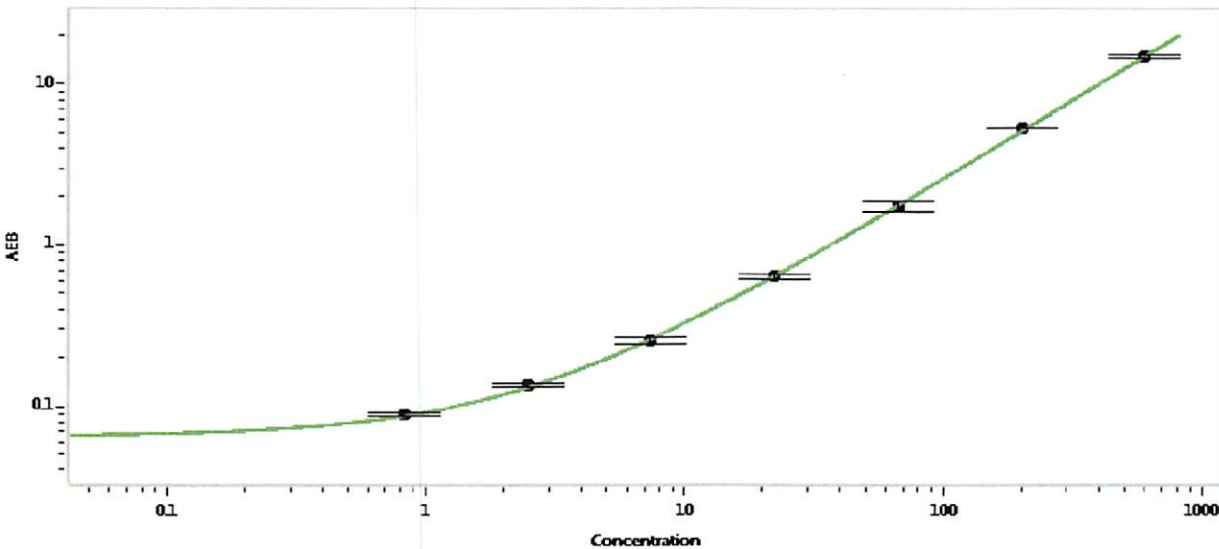
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102201	921405
Detector Stock	102202	921309
Calibrator Concentrate	102205	827108
Calibrator Diluent	102203	921207
Sample Diluent	102204	921208

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	12 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.823
C	2.47
D	7.41
E	22.2
F	66.7
G	200
H	600

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

Signature/Date