

Product Number:	102669
Lot Number:	502158
Expiration:	21-Nov-2020
Platform(s):	HD-1/HD-X



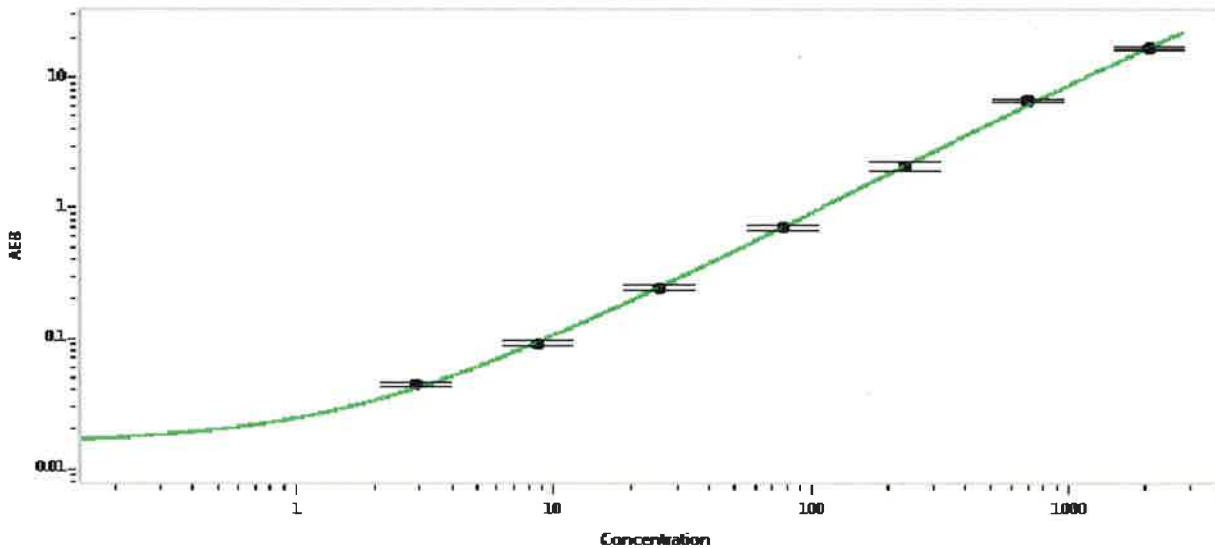
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102662	934410
Detector Stock	102663	934005
Calibrator Concentrate	102666	934501
Calibrator Diluent	102664	934022
Sample Diluent	102665	934006

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
Signature/Date
27 Jun 2020