

Product Number:	102648
Lot Number:	502014
Expiration:	30-Apr-2020
Platform(s):	HD-1/HD-X & SR-X

Component	Part Number	Lot Number
Bead Stock	102641	924609
Detector Stock	102642	924903
Detector Diluent	102734	924106
Calibrator Concentrate	102645	924602
Calibrator Diluent	102643	924007
Sample Diluent	102644	912024

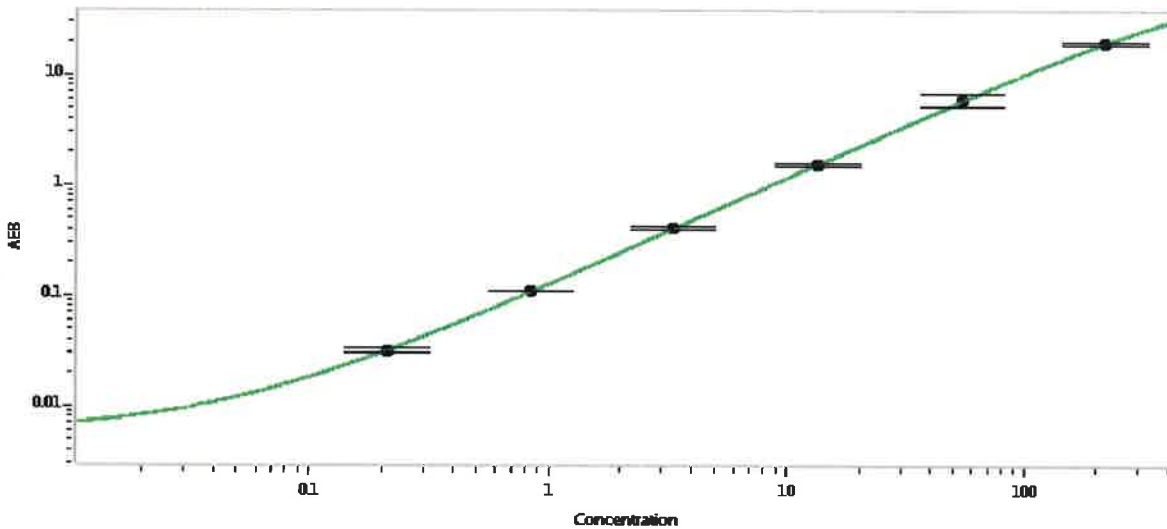


Data below represents results generated on the Simoa™ HD-1 Analyzer.

Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159

Calibrator Concentrate Target Value	3,000 pg/mL
-------------------------------------	-------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.21
C	0.84
D	3.36
E	13.4
F	53.8
G	215

Review/Approval: Patrizia Stadler Assoc. Director, QC P. Stadler 12 NOV 2019  
Name Title Signature/Date