

Product Number:	102794
Lot Number:	502294
Expiration:	11-Dec-2020
Platform(s):	HD-1/HD-X



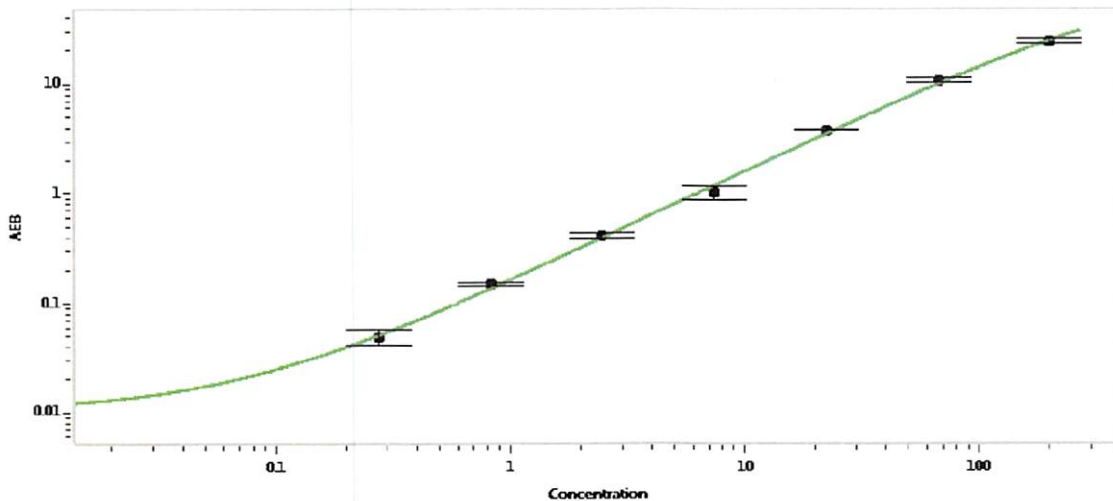
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102787	935111
Detector Stock	102788	935305
Calibrator Concentrate	102791	936007
Calibrator Diluent	102789	935328
Sample Diluent	102790	935320

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval

Patrizia Stadler  
Name

Assoc. Director, QC  
Title

*P. Stodler* 20 MAR 2020  
Signature/Date