

Product Number:	102517
Lot Number:	502337
Expiration:	19-Dec-2020
Platform(s):	HD-1/HD-X & SR-X



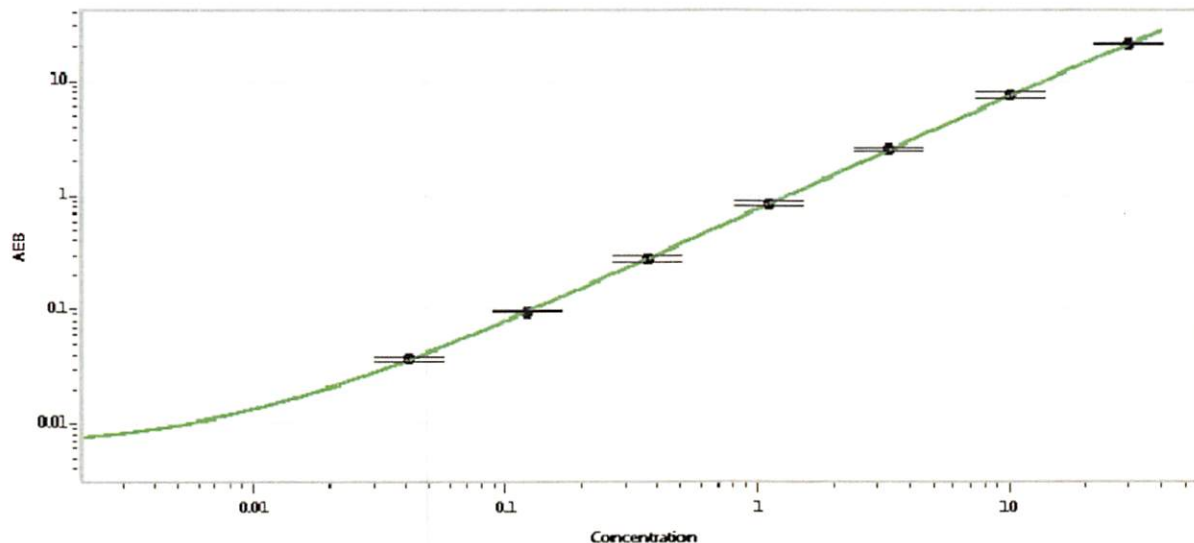
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102510	935327
Detector Stock	102511	935403
Calibrator Concentrate	102514	936010
Calibrator Diluent	102512	935331
Sample Diluent	102513	935324

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	600 pg/mL
-------------------------------------	-----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i> 21 Apr 2020
	Name	Title	Signature/Date