

Product Number:	103093
Lot Number:	502159
Expiration:	21-Nov-2020
Platform(s):	HD-1/ HD-X & SR-X



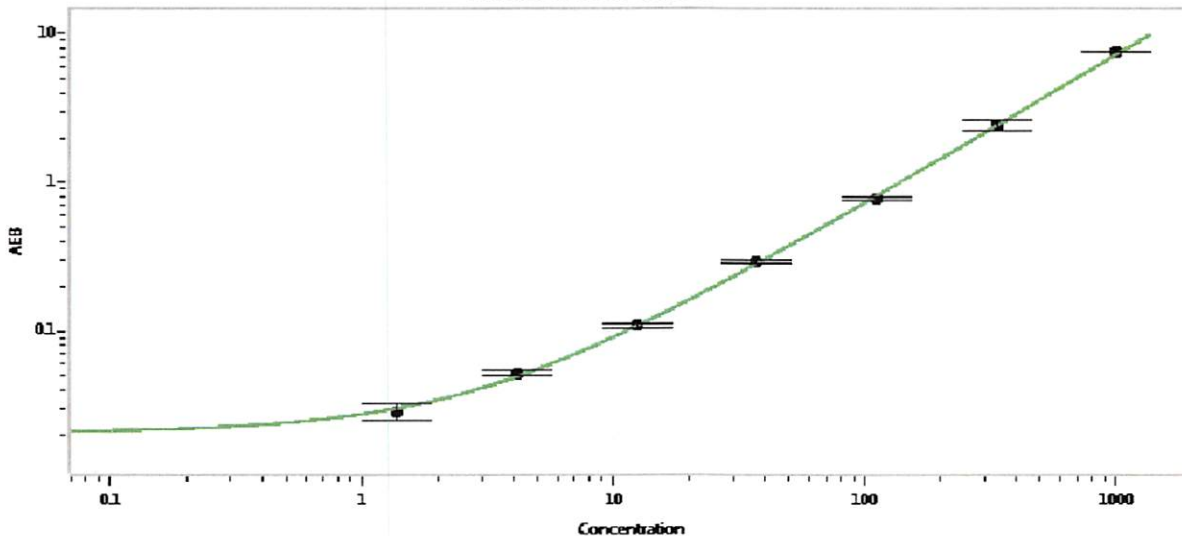
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102021	934609
Detector Stock	103043	935318
Calibrator Concentrate	103047	935308
Calibrator Diluent	103044	935307
Sample Diluent	103046	935319

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	1.37
C	4.12
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler 11/19/20</i>
	Name	Title	Signature/Date