

Product Number:	102336
Lot Number:	502142
Expiration:	1-Nov-2020
Platform(s):	HD-1/HD-X



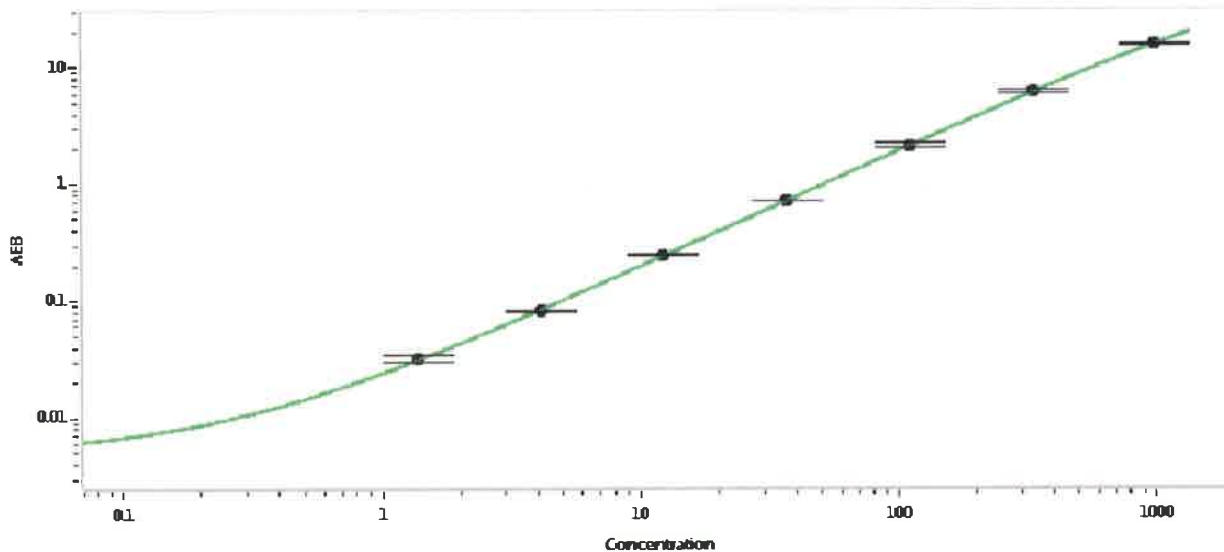
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102330	932914
Detector Stock	102331	932613
Calibrator Concentrate	102335	933104
Calibrator Diluent	102332	932615
Sample Diluent	102333	932617

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.37
C	4.11
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	16 Dec 2019
	Name	Title	Signature/Date	