

Product Number:	102491
Lot Number:	502143
Expiration:	27-May-2020
Platform(s):	HD-1/HD-X



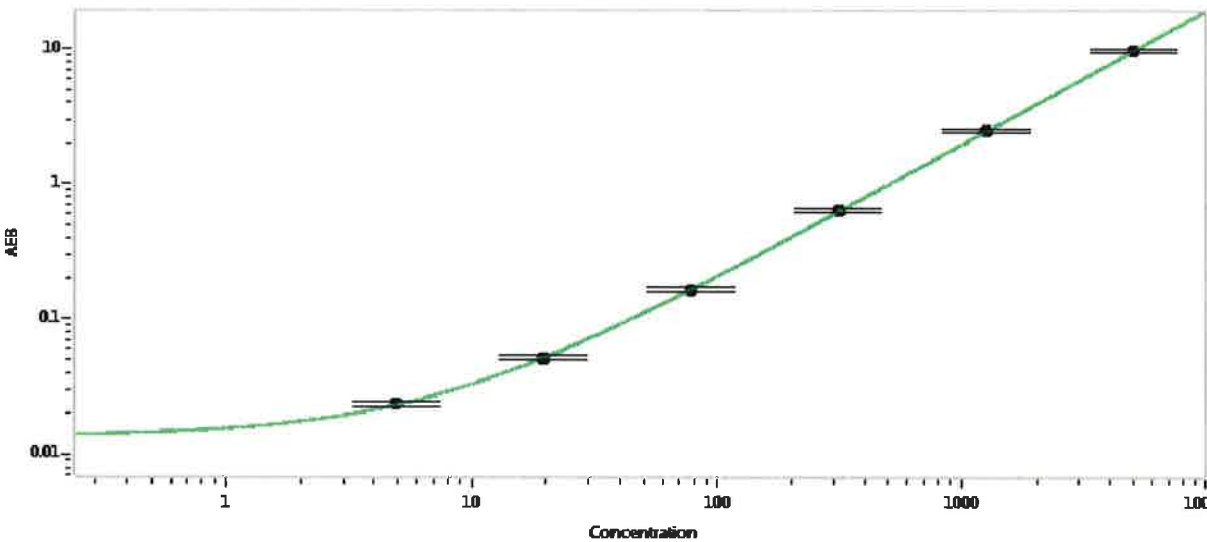
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102484	933805
Detector Stock	102485	933906
Detector Diluent	103084	933115
Calibrator Concentrate	102488	907004
Calibrator Diluent	102486	933101
Sample Diluent	102487	933102

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	4.88
C	19.5
D	78.1
E	313
F	1250
G	5000

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