

Product Number:	102073
Lot Number:	501967
Expiration:	13-Nov-2019
Platform(s):	HD-1/HD-X



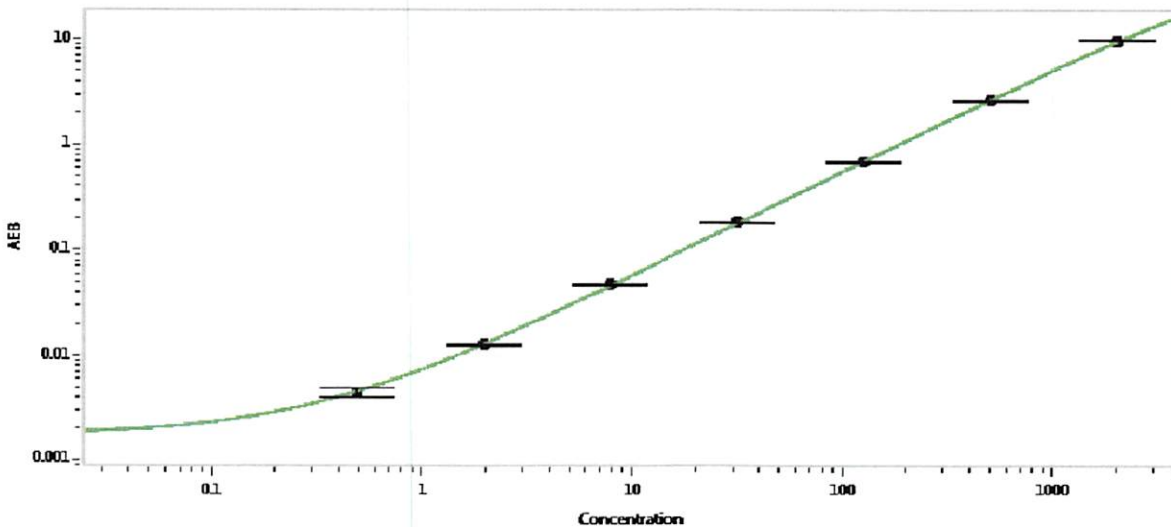
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102065	835109
Detector Stock	102066	902501
Calibrator Concentrate	102069	834103
Calibrator Diluent	102067	831713
Sample Diluent	102068	831714

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

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.488
C	1.95
D	7.81
E	31.3
F	125
G	500
H	2000

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P Stadler</i> 02 Aug 2019
	Name	Title	Signature/Date