

Product Number:	102570
Lot Number:	502064
Expiration:	1-Aug-2020
Platform(s):	HD-1/HD-X



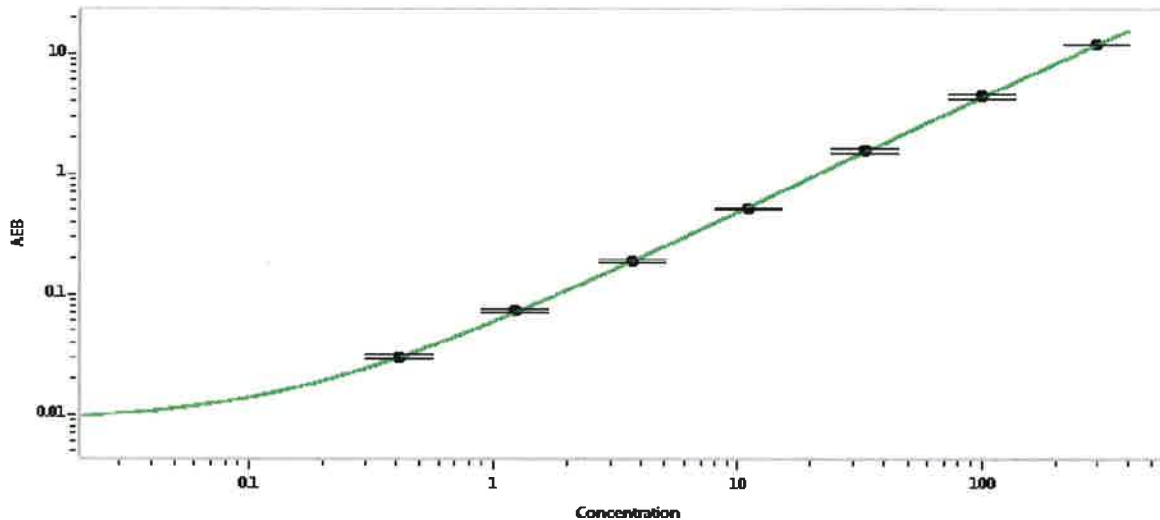
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102563	928101
Detector Stock	102564	927711
Detector Diluent	102590	928208
Calibrator Concentrate	102567	829005
Calibrator Diluent	102565	927713
Sample Diluent	102566	927714

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	6 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.412
C	1.23
D	3.70
E	11.1
F	33.3
G	100
H	300

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	18 Oct 2019
	Name	Title	Signature/Date	