

Product Number:	101984
Lot Number:	502007
Expiration:	25-Jan-2020
Platform(s):	HD-1/HD-X



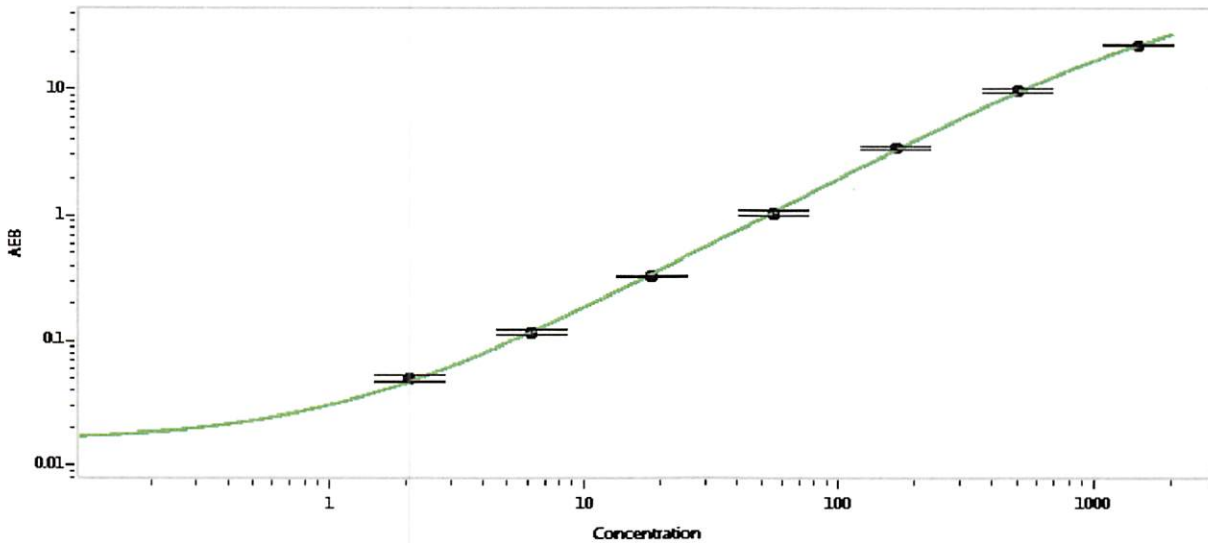
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	101910	902504
Detector Stock	101911	909115
Calibrator Concentrate	101983	910010
Calibrator Diluent	101908	902603
Sample Diluent	101909	908505

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	30 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	2.06
C	6.2
D	18.5
E	56
F	167
G	500
H	1500

Review/Approval	Patrizia Stadler	QC Assoc. Director	<i>P. Stadler</i>	06 Sep 2019
	Name	Title	Signature/Date	