

Product Number:	101984
Lot Number:	502074
Expiration:	24-Sep-2020
Platform(s):	HD-1/HD-X



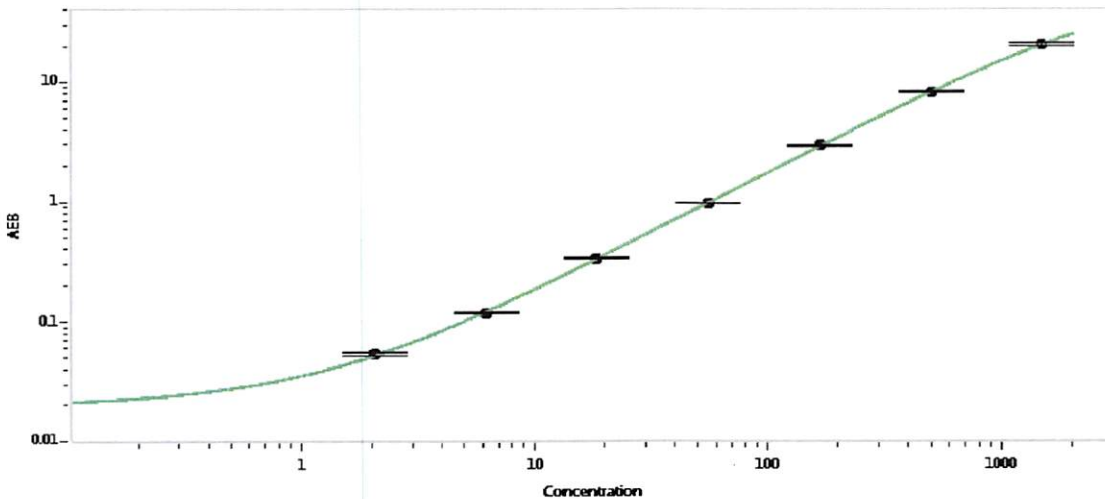
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	101910	928103
Detector Stock	101911	928010
Calibrator Concentrate	101983	910010
Calibrator Diluent	101908	928203
Sample Diluent	101909	928011

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	6.17
D	18.5
E	55.6
F	167
G	500
H	1500

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>p.stadler</i>	17 Oct 2019
	Name	Title	Signature/Date	