

Product Number:	102318
Lot Number:	501802
Expiration:	3-Dec-2019
Platform(s):	HD-1



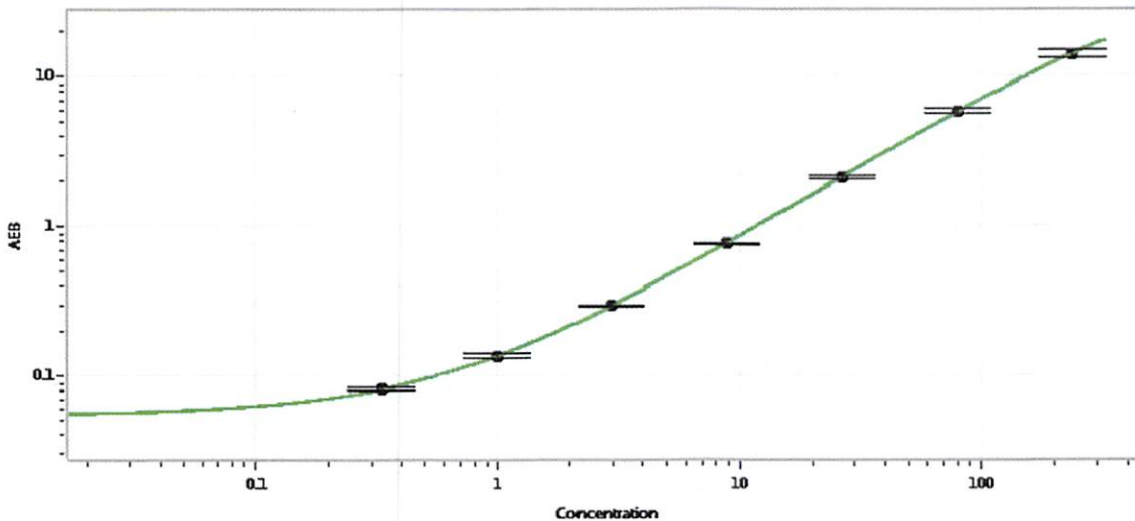
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102309	912001
Detector Stock	102310	908520
Calibrator Concentrate	102314	908607
Calibrator Diluent	102311	833714
Sample Diluent	102312	833713

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	4,800 pg/mL
-------------------------------------	-------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.329
C	0.988
D	2.96
E	8.89
F	26.7
G	80.0
H	240

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	02/17/2019
	Name	Title	Signature/Date	