

Product Number:	102318
Lot Number:	501654
Expiration:	20-Nov-2019
Platform(s):	HD-1



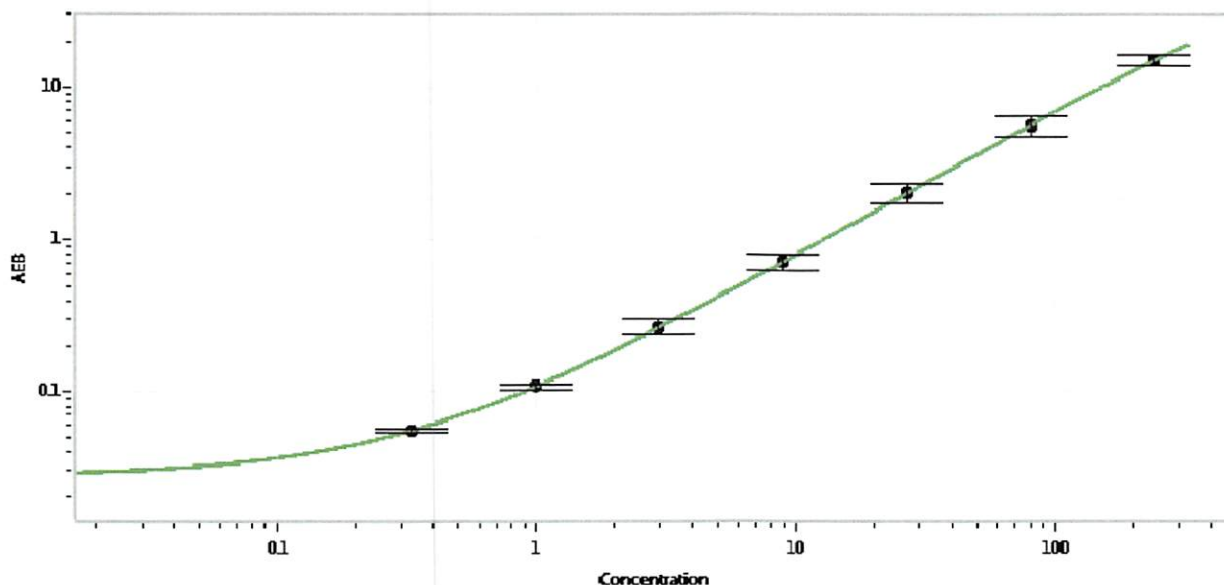
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102309	833804
Detector Stock	102310	834104
Calibrator Concentrate	102314	818414
Calibrator Diluent	102311	832415
Sample Diluent	102312	832416

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> 14 Feb 2019
	Name	Title	Signature/Date