

<b>Product Number:</b>	102336
<b>Lot Number:</b>	501726
<b>Expiration:</b>	21-Dec-2019
<b>Platform(s):</b>	HD-1

Component	Part Number	Lot Number
Bead Stock	102330	908105
Detector Stock	102331	907303
Calibrator Concentrate	102335	826303
Calibrator Diluent	102332	835507
Sample Diluent	102333	907011

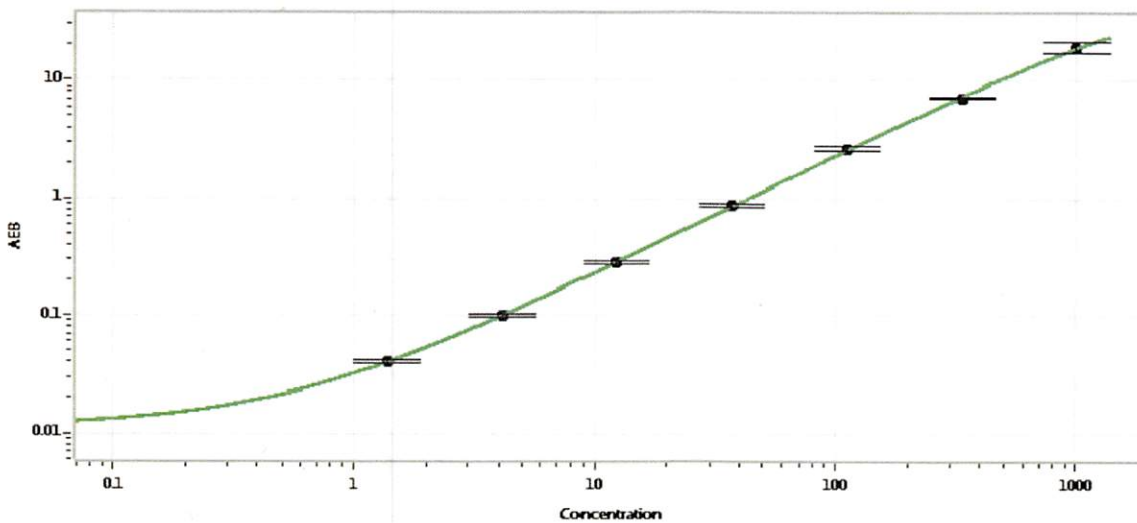
Generic Components - Included With Kit But Not Lot Number Specific	
Discovery Bead Diluent	101954
Discovery Detector Diluent	101955
Discovery SBG Diluent	101956
Discovery SBG Stock	103297
RGP Reagent	103159



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

<b>Calibrator Concentrate Target Value</b>	20 ng/mL
--	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.37
C	4.12
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval

Patrizia Stadler  
Name

QC Assoc. Director  
Title

*P. Stadler* 28 Mar 2019  
Signature/Date