

Product Number:	102635
Lot Number:	502228
Expiration:	9-Aug-2020
Platform(s):	HD-1/ HD-X

Component	Part Number	Lot Number
Bead Stock	102628	002811
Detector Stock	102629	003604
Calibrator Concentrate	102632	822101
Calibrator Diluent	102630	002216
Sample Diluent	102631	002804

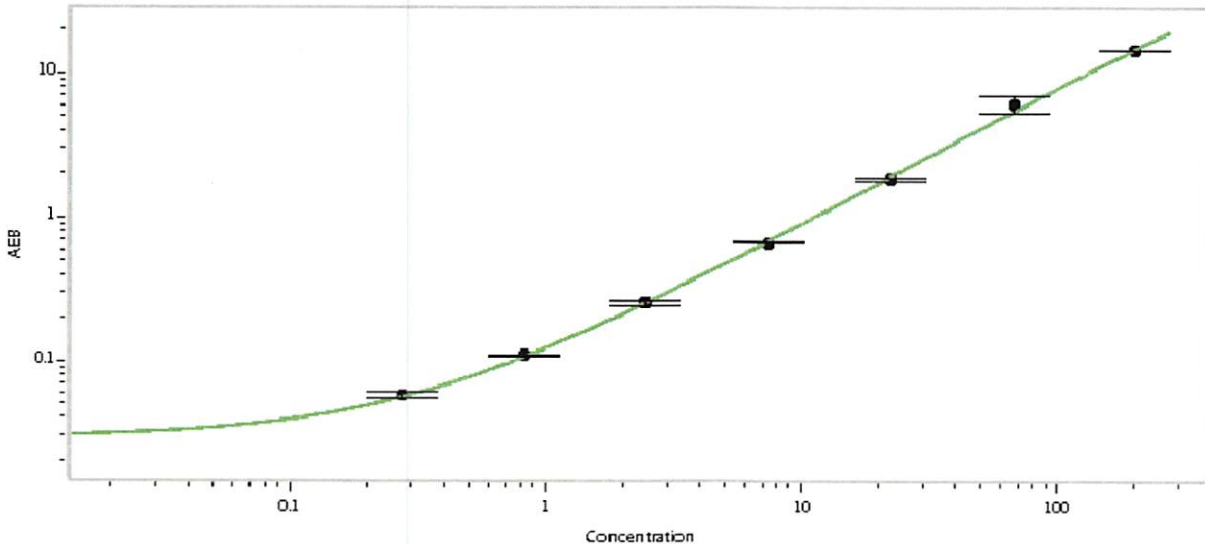
Generic Components - Included With Kit But Not Lot Number Specific		
Discovery Bead Diluent		101954
Discovery Detector Diluent		101955
Discovery SBG Diluent		101956
Discovery SBG Stock		103297
RGP Reagent		103159



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

Calibrator Concentrate Target Value	4 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler 06 Feb 2020
Signature/Date