

Product Number:	102039
Lot Number:	501656
Expiration:	24-Oct-2019
Platform(s):	HD-1 & SR-X

Component	Part Number	Lot Number
Bead Stock	102028	832315
Detector Stock	102029	833109
Calibrator Concentrate	102032	MF-042618-03
Calibrator Diluent	102030	829712
Sample Diluent	102031	831816

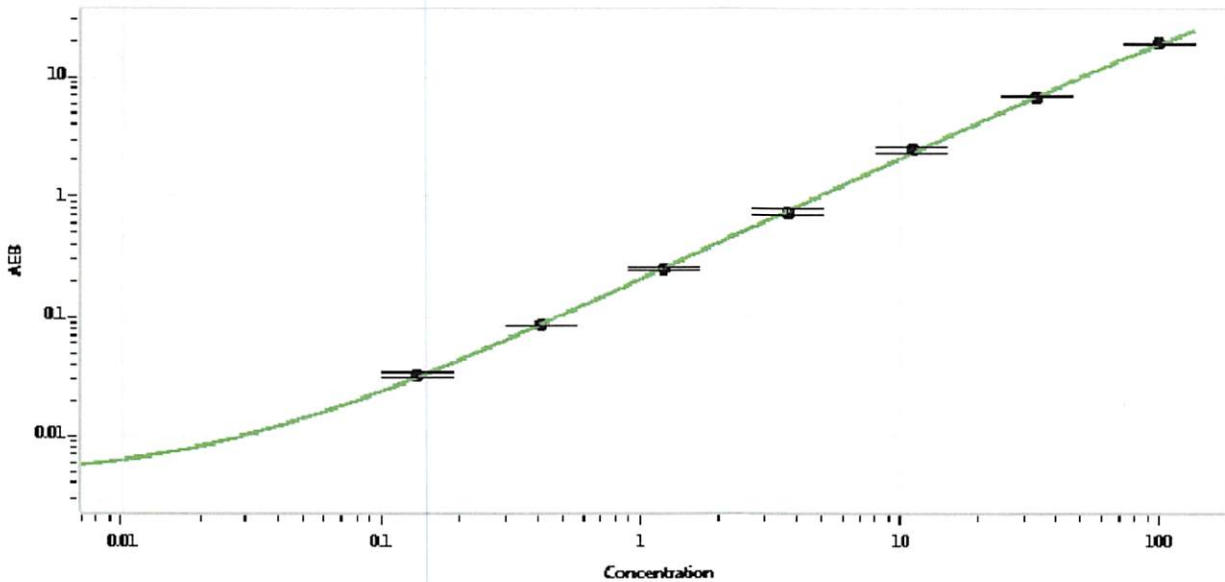
Generic Components - Included With Kit But Not Lot Number Specific		
Discovery Bead Diluent		101954
Discovery Detector Diluent		101955
Discovery SBG Diluent		101956
Discovery SBG Stock		103297
RGP Reagent		103159



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

Calibrator Concentrate Target Value	2 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.137
C	0.412
D	1.23
E	3.70
F	11.1
G	33.3
H	100

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler 12 Feb 2019
Signature/Date