

Product Number:	102394
Lot Number:	501976
Expiration:	5-Jan-2020
Platform(s):	HD-1/HD-X

Component	Part Number	Lot Number
Bead Stock	102387	921804
Detector Stock	102388	921805
Calibrator Concentrate	102391	MF-010518-06
Calibrator Diluent	102389	919311
Sample Diluent	102390	912021

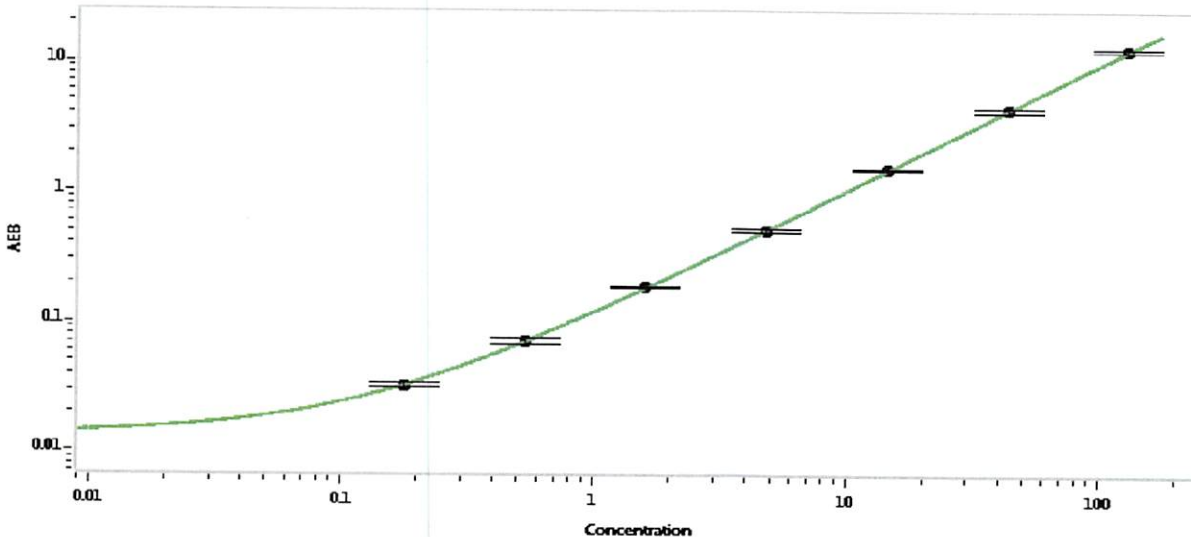


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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	4000 pg/mL
-------------------------------------	------------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.178
C	0.535
D	1.60
E	4.81
F	14.4
G	43.3
H	130

Review/Approval

Patrizia Stadler  
Name

Assoc. Director, QC  
Title

*P. Stadler* 09 Aug 2019  
Signature/Date