

Product Number:	102794
Lot Number:	501735
Expiration:	19-Jan-2020
Platform(s):	HD-1

Component	Part Number	Lot Number
Bead Stock	102787	903204
Detector Stock	102788	903108
Calibrator Concentrate	102791	MF-021018-04
Calibrator Diluent	102789	901914
Sample Diluent	102790	903203

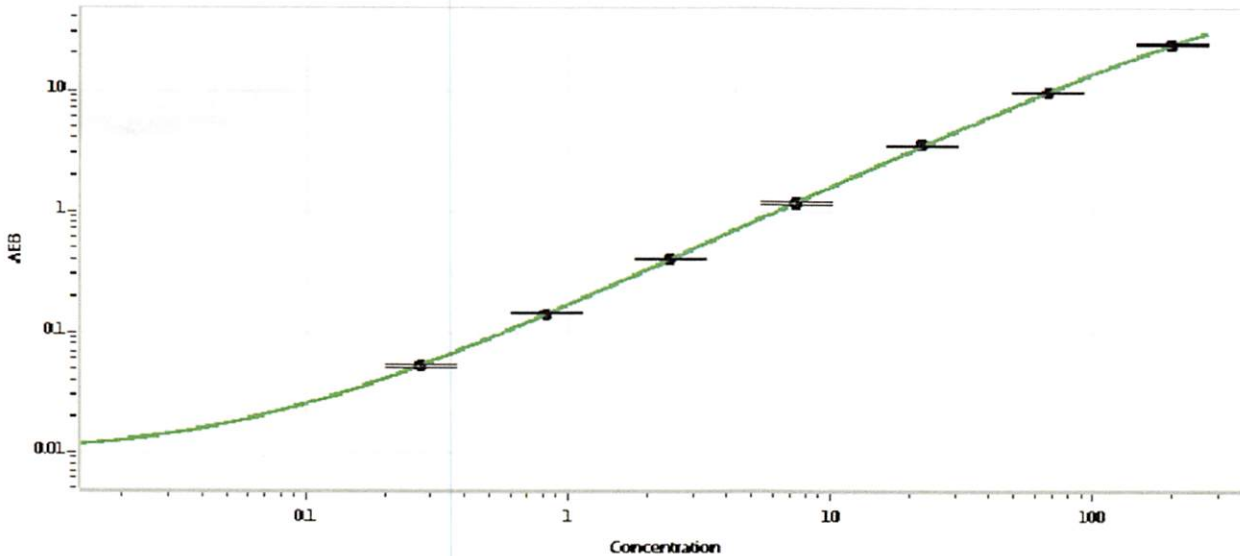


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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	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 22 MAR 2019
Signature/Date