

Product Number:	102064
Lot Number:	502448
Expiration:	3-Dec-2020
Platform(s):	HD-1/HD-X

Component	Part Number	Lot Number
Bead Stock	102056	018204
Detector Stock	102057	018102
Calibrator Concentrate	102060	833709
Calibrator Diluent	102058	018207
Sample Diluent	102059	018206

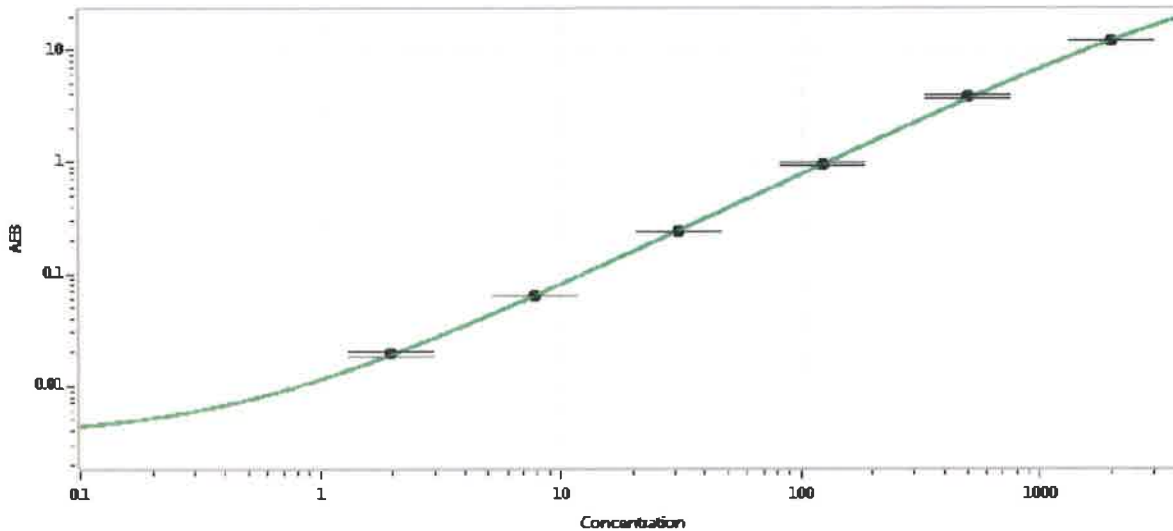


Data below represents results generated on the Simoa™ HD-1 Analyzer.

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	40 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.95
C	7.81
D	31.3
E	125
F	500
G	2000

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P Stadler 17 Jul 2020
Signature/Date