

Product Number:	102700
Lot Number:	502197
Expiration:	4-Jun-2020
Platform(s):	HD-1/HD-X



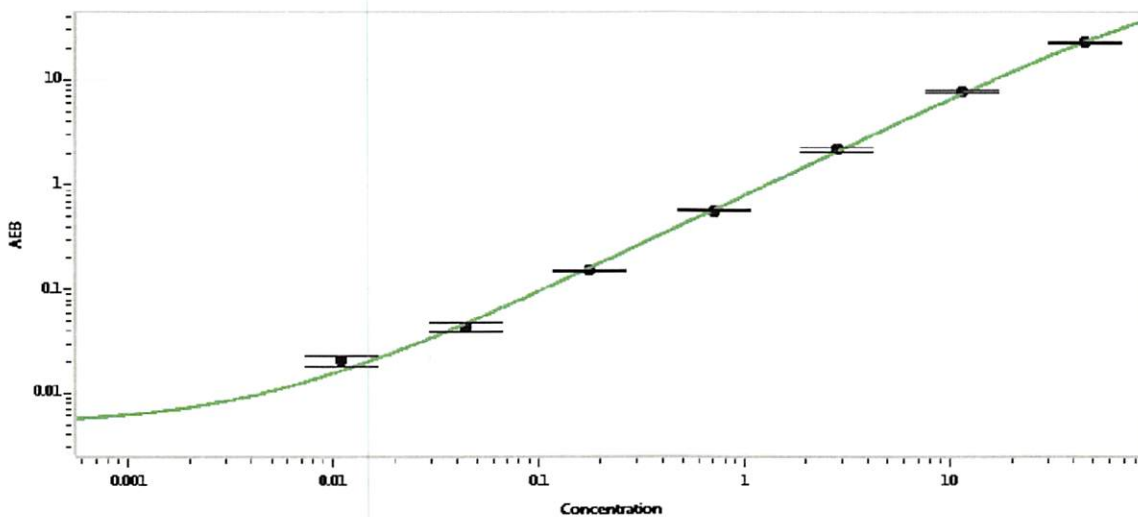
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102693	001306
Detector Stock	102694	001310
Calibrator Concentrate	102697	MF-060418-01
Calibrator Diluent	102695	933811
Sample Diluent	102696	933812

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	1000 pg/mL
-------------------------------------	------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.00
B	0.011
C	0.044
D	0.176
E	0.703
F	2.81
G	11.3
H	45.0

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler 16 Jan 2020
Signature/Date