

Product Number:	101419
Lot Number:	502193
Expiration:	13-Oct-2020
Platform(s):	HD-1/HD-X



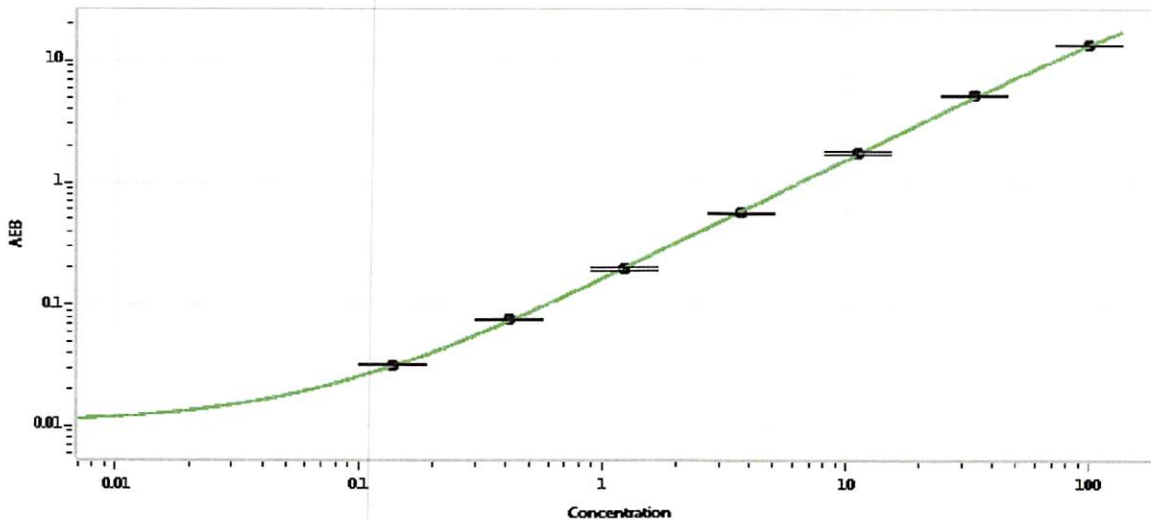
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	101899	001307
Detector Stock	101900	001311
Calibrator Concentrate	101420	828601
Calibrator Diluent	101418	932924
Sample Diluent	101416	932624

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	2 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.137
C	0.412
D	1.24
E	3.70
F	11.1
G	33.3
H	100

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler 17 Jan 2020
Signature/Date