

Product Number:	102318
Lot Number:	501743
Expiration:	11-Feb-2020
Platform(s):	HD-1



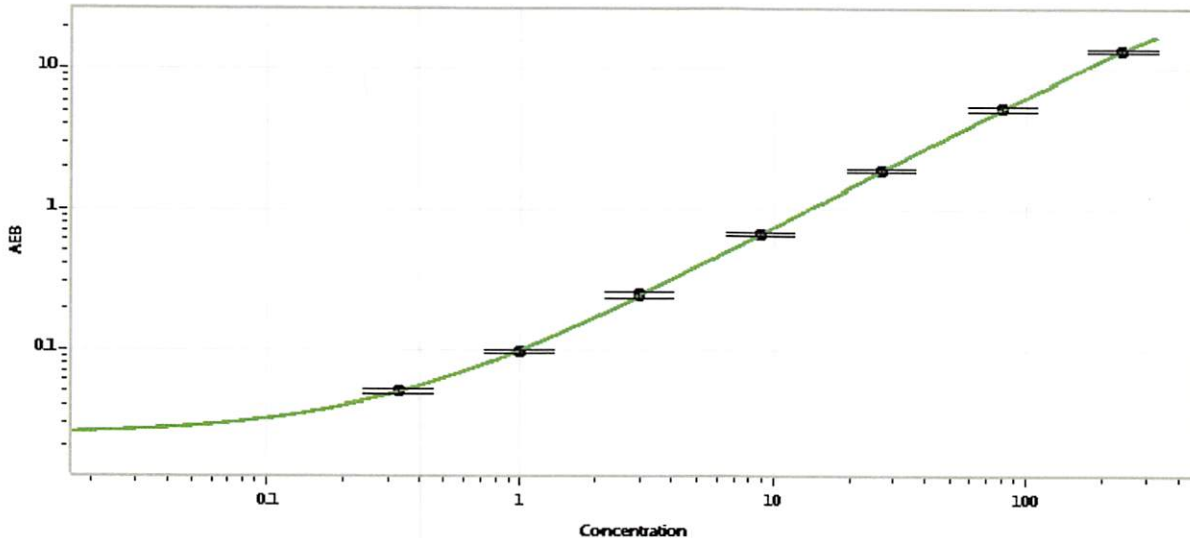
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102309	908506
Detector Stock	102310	908520
Calibrator Concentrate	102314	908607
Calibrator Diluent	102311	907703
Sample Diluent	102312	907704

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	4,800 pg/mL
-------------------------------------	-------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.329
C	0.988
D	2.96
E	8.89
F	26.7
G	80.0
H	240

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler 29 Mar 2017
Signature/Date