

Product Number:	102318
Lot Number:	501891
Expiration:	26-Mar-2020
Platform(s):	HD-1/HD-x



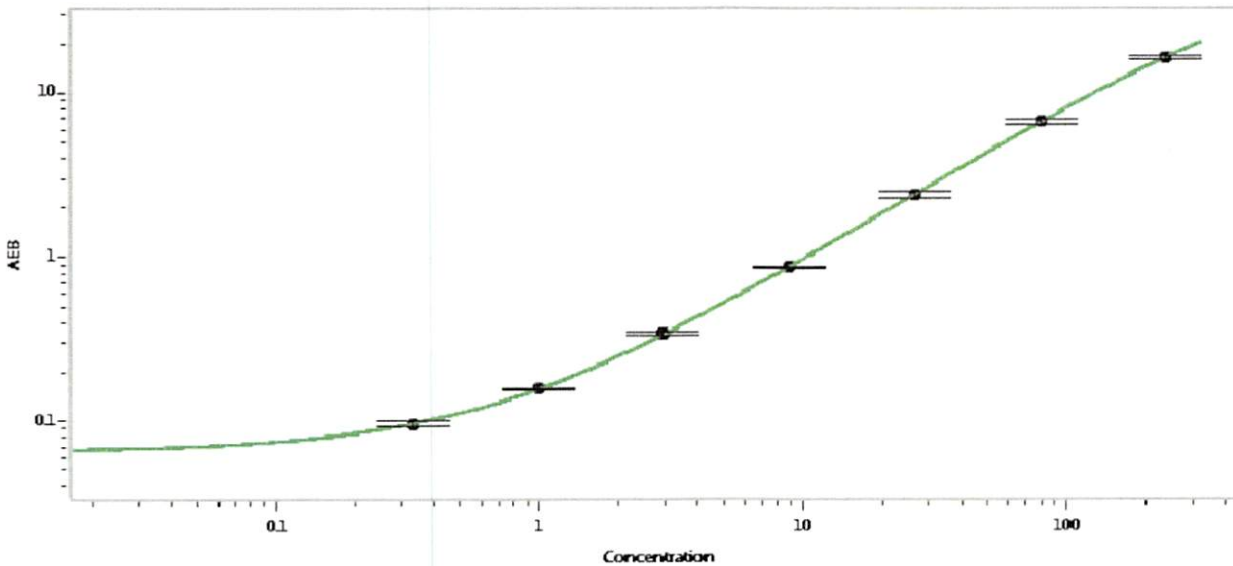
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102309	918211
Detector Stock	102310	908520
Calibrator Concentrate	102314	908607
Calibrator Diluent	102311	917904
Sample Diluent	102312	917905

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 03 JUL 2019
Signature/Date