

Product Number:	102343
Lot Number:	502030
Expiration:	11-Mar-2020
Platform(s):	HD-1/HD-X



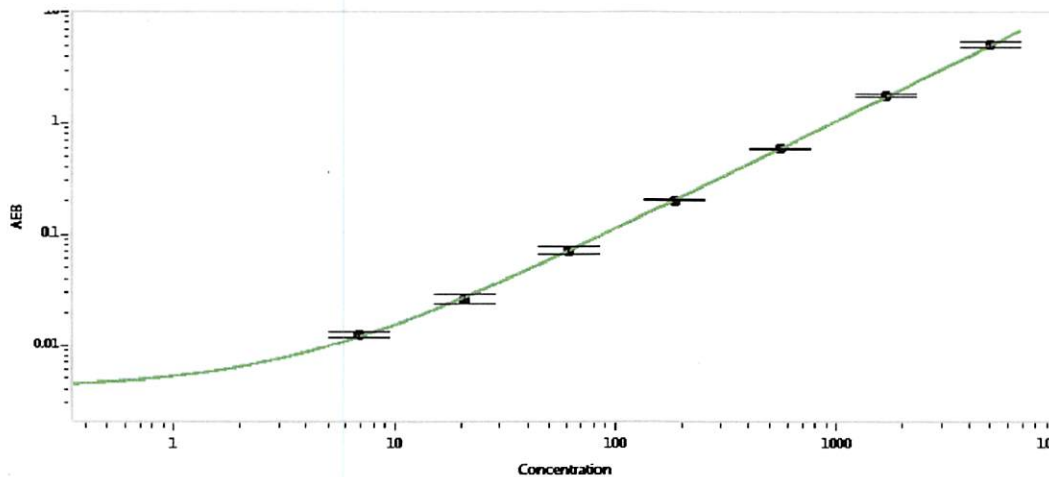
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102337	907707
Detector Stock	102338	925902
Calibrator Concentrate	102342	831815
Calibrator Diluent	102339	908919
Sample Diluent	102340	907008

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	100 ng/mL
-------------------------------------	-----------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.00
B	6.86
C	20.6
D	61.7
E	185
F	556
G	1667
H	5000

Review/Approval

Patrizia Stadler  
Name

Assoc. Director, QC  
Title

*P. Stadler 25 Sep 2019*  
Signature/Date