

Product Number:	102336
Lot Number:	502046
Expiration:	30-Mar-2020
Platform(s):	HD-1/HD-X



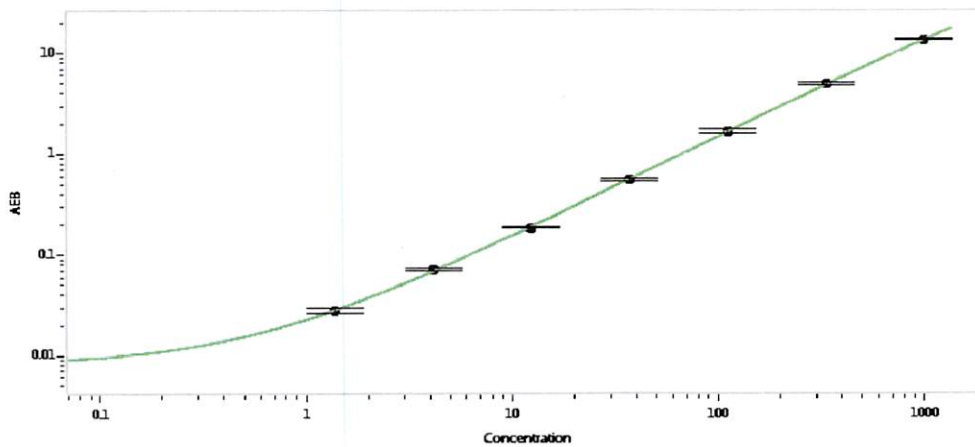
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102330	926607
Detector Stock	102331	926606
Calibrator Concentrate	102335	926406
Calibrator Diluent	102332	908922
Sample Diluent	102333	925304

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	20 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.37
C	4.11
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler 27 Sep 2019
Signature/Date