

Product Number:	102336
Lot Number:	501618
Expiration:	1-Oct-2019
Platform(s):	HD-1



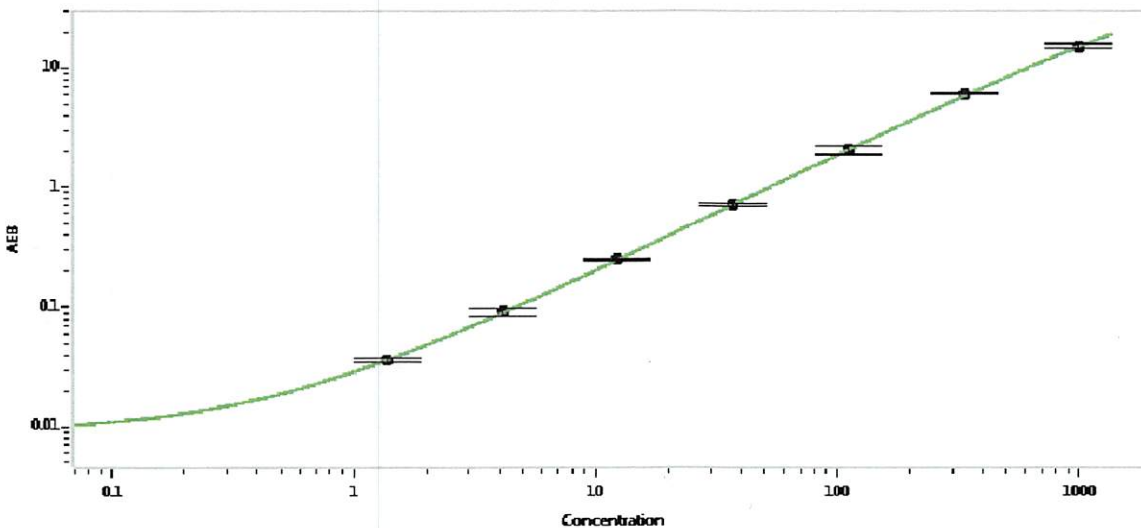
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102330	902907
Detector Stock	102331	902801
Calibrator Concentrate	102335	826303
Calibrator Diluent	102332	827416
Sample Diluent	102333	902208

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 QC Assoc. Director
Name Title

P. Stodler 01 Feb 2019
Signature/Date