

Product Number:	102184
Lot Number:	501968
Expiration:	30-Apr-2020
Platform(s):	HD-1/ HD-X



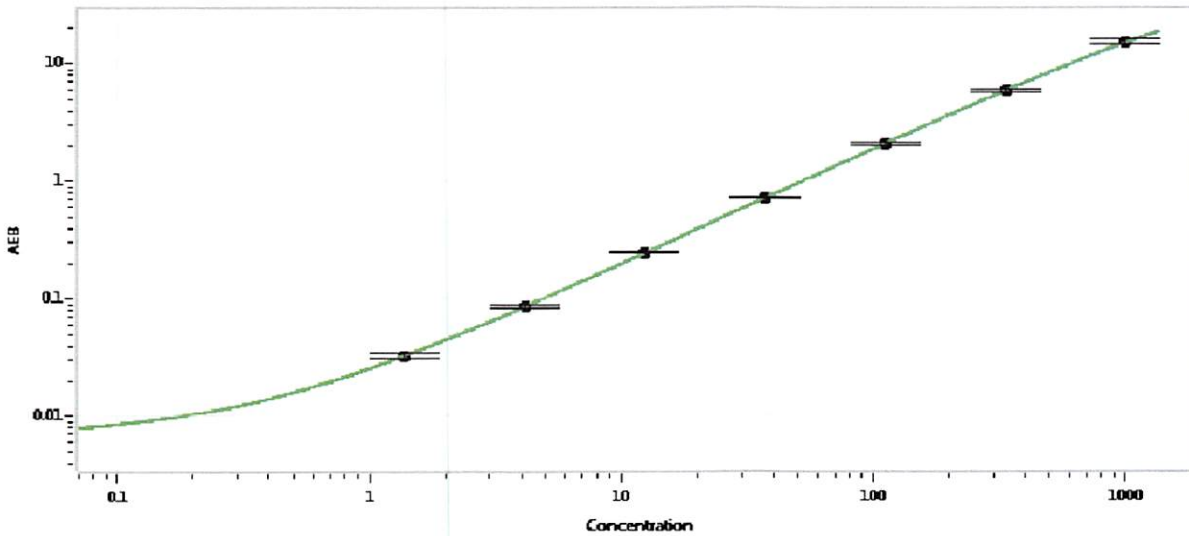
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102176	921010
Detector Stock	102177	921307
Calibrator Concentrate	102180	830509
Calibrator Diluent	102178	919310
Sample Diluent	102179	912020

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	40 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	1.37
C	4.12
D	12.3
E	37.0
F	111
G	333
H	1000

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler 08 Aug 2019
Signature/Date