

Product Number:	102713
Lot Number:	501914
Expiration:	30-Apr-2020
Platform(s):	HD-1/HD-X



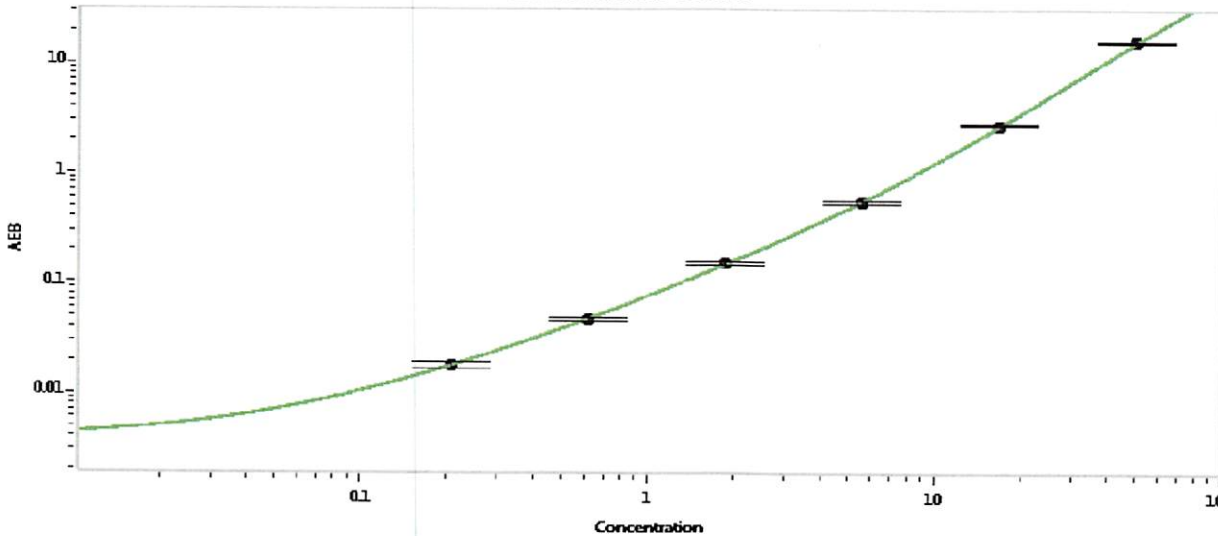
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102706	919608
Detector Stock	102707	921707
Calibrator Concentrate	102710	919307
Calibrator Diluent	102708	919009
Sample Diluent	102709	912019

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	2 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.206
C	0.617
D	1.85
E	5.56
F	16.7
G	50.0

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler 07 Aug 2019
Signature/Date