

Product Number:	102184
Lot Number:	501624
Expiration:	24-Aug-2019
Platform(s):	HD-1



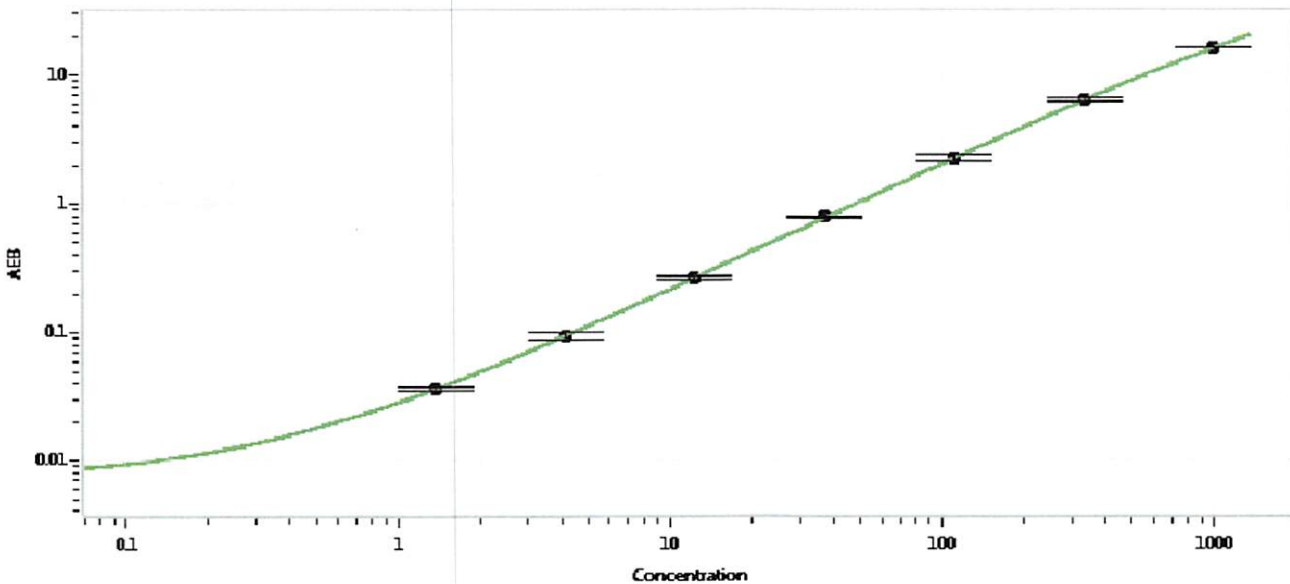
Data below represents results generated on the Simoa™ HD-1

Component	Part Number	Lot Number
Bead Stock	102176	827508
Detector Stock	102177	828311
Calibrator Concentrate	102180	830509
Calibrator Diluent	102178	823614
Sample Diluent	102179	826209

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

QC Assoc. Director
Title

P. Stadler 29 Jan 2019
Signature/Date