

Product Number:	102073
Lot Number:	501616
Expiration:	2-Nov-2019
Platform(s):	HD-1



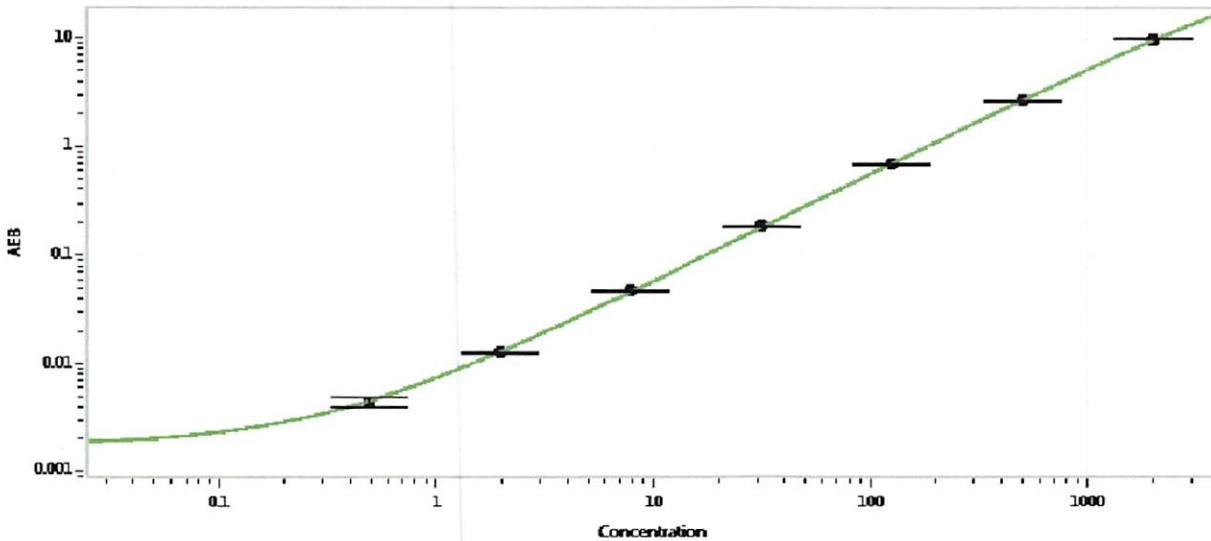
Data below represents results generated on the Simoa™ HD-1

Component	Part Number	Lot Number
Bead Stock	102065	835109
Detector Stock	102066	902501
Calibrator Concentrate	102069	834103
Calibrator Diluent	102067	831713
Sample Diluent	102068	831714

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	0.488
C	1.95
D	7.81
E	31.3
F	125
G	500
H	2000

Review/Approval

Patrizia Stadler
Name

QC Assoc. Director
Title

P. Stadler 12Feb2019
Signature/Date