

Product Number:	102669
Lot Number:	501615
Expiration:	3-Dec-2019
Platform(s):	HD-1



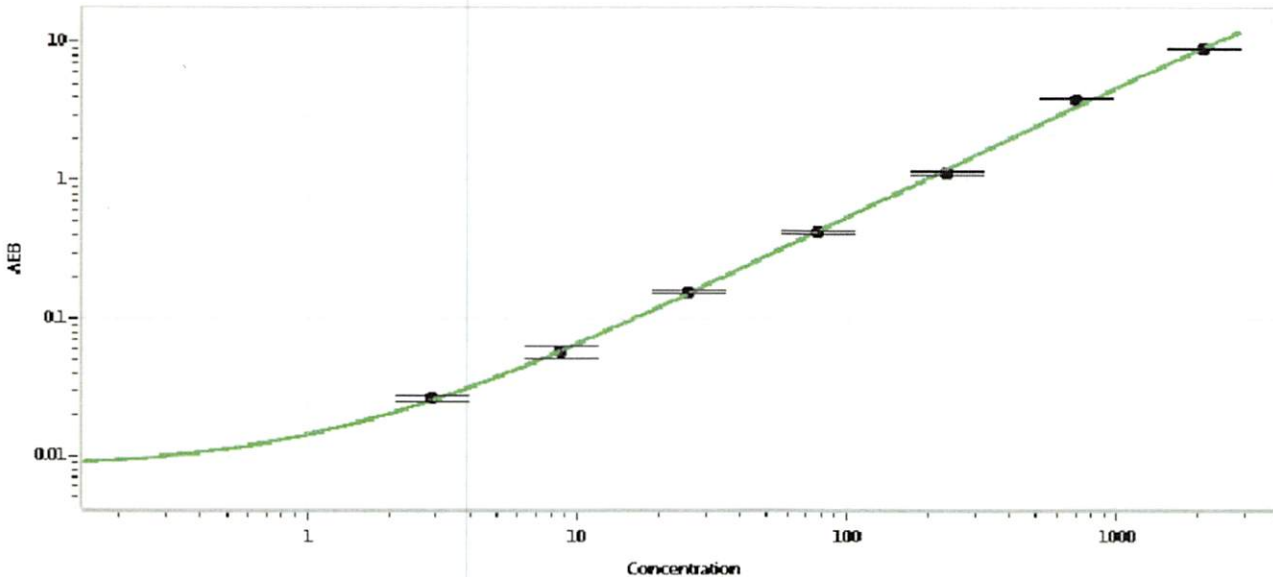
Data below represents results generated on the Simoa™ HD-1

Component	Part Number	Lot Number
Bead Stock	102662	902405
Detector Stock	102663	902505
Calibrator Concentrate	102666	MF-020718-03
Calibrator Diluent	102664	901911
Sample Diluent	102665	902506

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	60 ng/mL
-------------------------------------	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.88
C	8.64
D	25.9
E	77.8
F	233
G	700
H	2100

Review/Approval	Patrizia Stadler	QC Assoc. Director	<i>P. Stadler</i>	01Feb2019
	Name	Title	Signature/Date	