

Product Number:	101832
Lot Number:	501827
Expiration:	14-Mar-2020
Platform(s):	HD-1



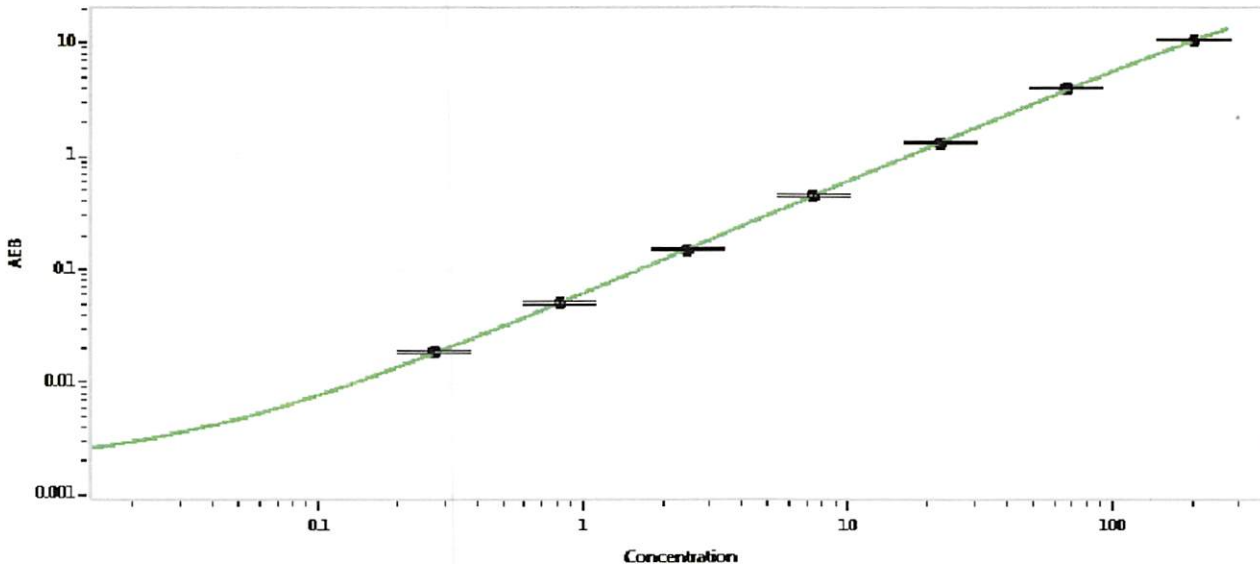
Data below represents results generated on the Simoa™ HD-1

Component	Part Number	Lot Number
Bead Stock	101927	912204
Detector Stock	101928	912806
Calibrator Concentrate	101827	MF-041618-09
Calibrator Diluent	101945	910311
Sample Diluent	101946	907317

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	4000 pg/mL
--	------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval	Patrizia Stadler	QC Assoc. Director	<i>p stadler</i>	15/04/2019
	Name	Title	Signature/Date	