

Product Number:	102475
Lot Number:	501800
Expiration:	14-Mar-2020
Platform(s):	HD-1



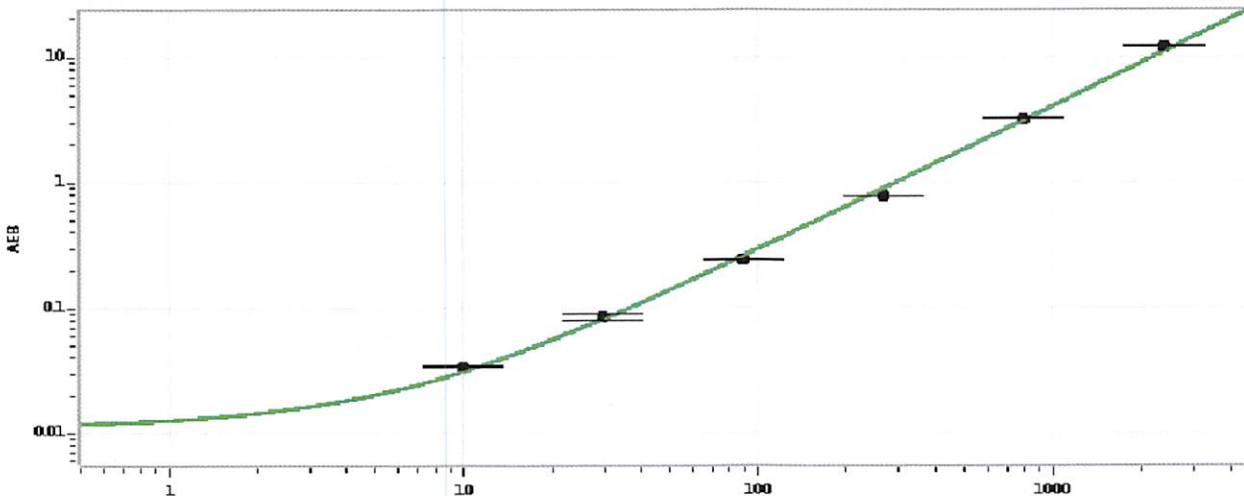
Data below represents results generated on the Simoa™ HD-1

Component	Part Number	Lot Number
Bead Stock	102468	911701
Detector Stock	102469	911702
Calibrator Concentrate	102472	MF-062018-06
Calibrator Diluent	102470	910309
Sample Diluent	102471	907315

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	9.88
C	29.6
D	88.9
E	267
F	800
G	2400

Review/Approval	Patrizia Stadler	QC Assoc. Director	<i>P. Stadler</i>	02 MAY 2019
	Name	Title	Signature/Date	