

Product Number:	102929
Lot Number:	502022
Expiration:	1-Aug-2020
Platform(s):	HD-1/HD-X



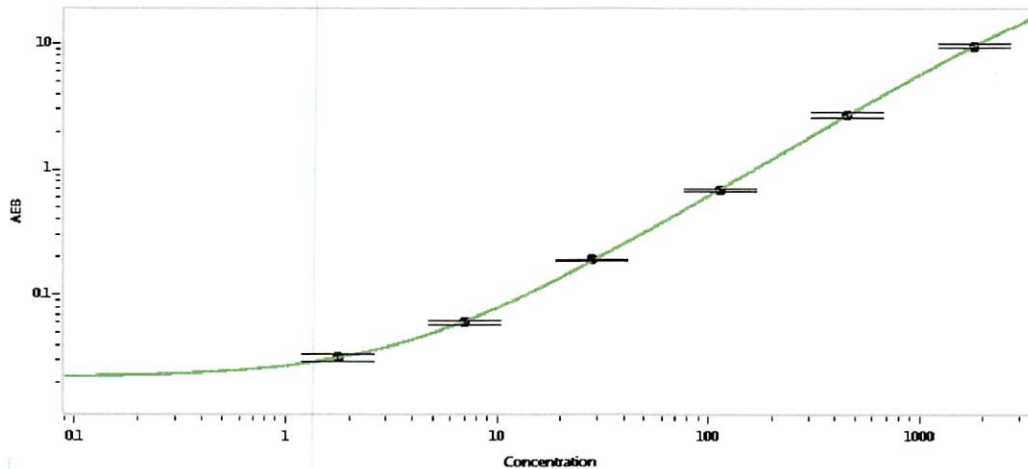
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102922	926312
Detector Stock	102923	926311
Calibrator Concentrate	102926	926313
Calibrator Diluent	102924	926310
Sample Diluent	102925	926210

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

<b>Calibrator Concentrate Target Value</b>	36 ng/mL
--	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	1.76
C	7.03
D	28.1
E	113
F	450
G	180

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>p. Stadler</i> 27 Sep 2019
	Name	Title	Signature/Date