

Product Number:	102039
Lot Number:	502622
Expiration:	16-Oct-2021
Platform(s):	HD-1/HD-X & SR-X



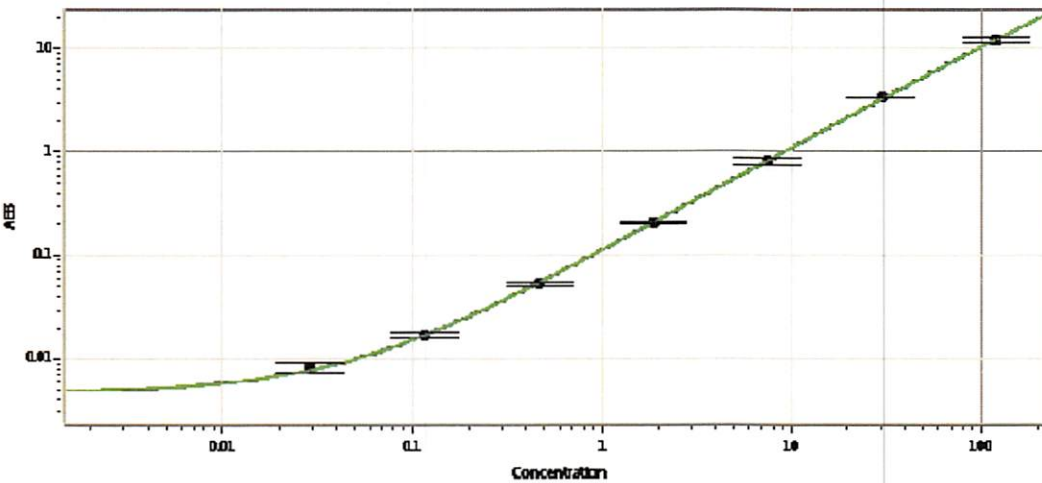
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102028	032207
Detector Stock	102029	032407
Calibrator Concentrate	102032	032408
Calibrator Diluent	102030	031412
Sample Diluent	103402	030405

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	2 ng/mL
-------------------------------------	---------

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.029
C	0.117
D	0.469
E	1.88
F	7.50
G	30.0
H	120

Review/Approval

Patrizia Stadler

Associate Director, QC

P. Stadler 25 Nov 2016

Name

Title

Signature/Date