

Product Number:	102233
Lot Number:	502265
Expiration:	24-Aug-2020
Platform(s):	HD-1/HD-X, SRX



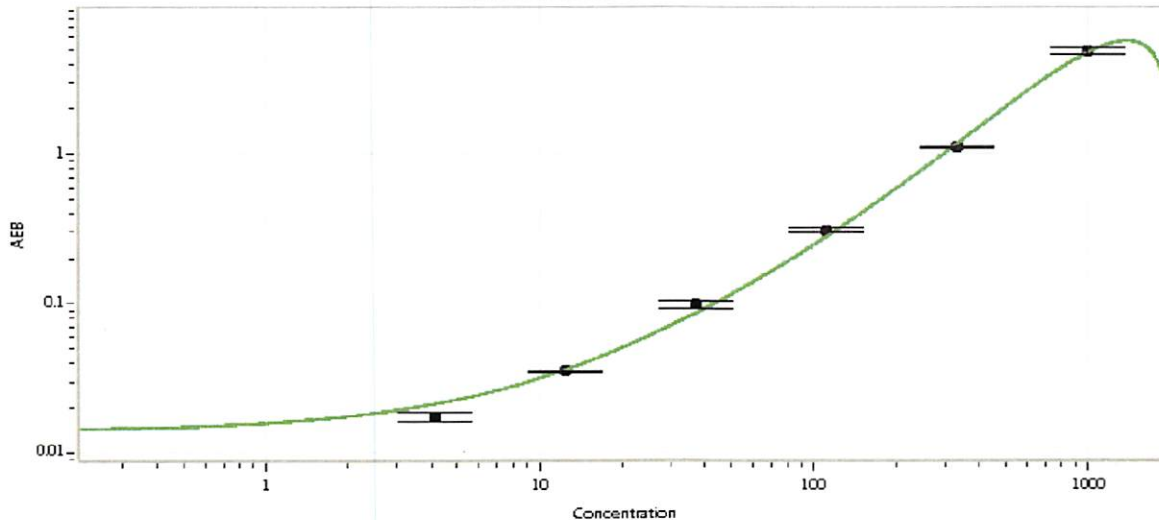
Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Stock	102223	005510
Detector Stock	102224	005511
Calibrator Concentrate	102229	835002
Calibrator Diluent	102225	005907
Sample Diluent	102226	005604

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	20 ng/mL
--	----------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	4.12
C	12.3
D	37.0
E	111
F	333
G	1000

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler 04/08/20

Name

Title

Signature/Date